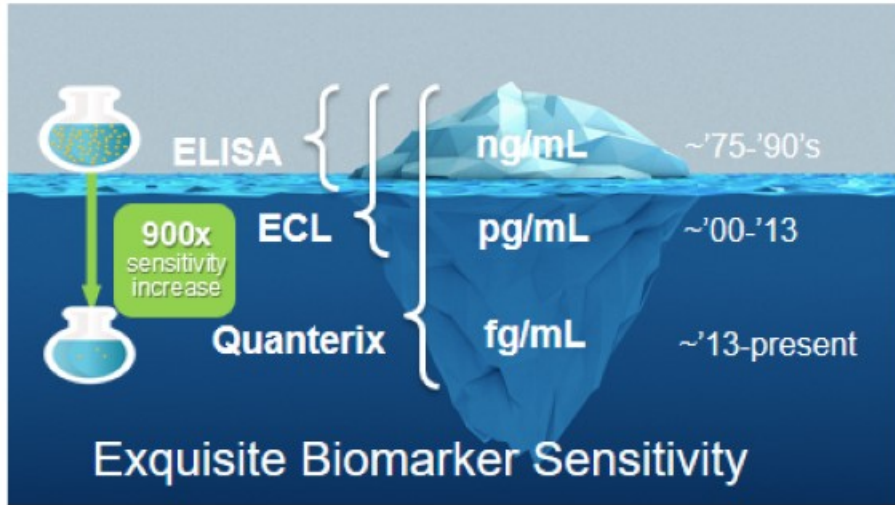


Q1 2019 Earnings Call • May 9, 2019

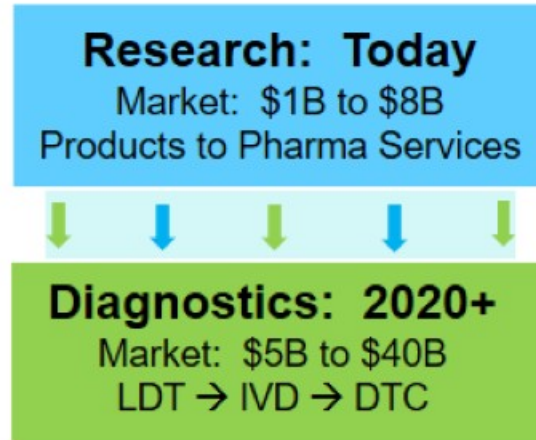
This presentation contains “forward-looking” statements that are based on our beliefs and assumptions and on information available to us as of the date of this presentation. Forward-looking statements include all statements that are not historical facts. We may not actually achieve the plans, intentions or expectations disclosed in our forward-looking statements. Forward-looking statements involve known and unknown risks, uncertainties, assumptions and other factors that may cause our actual results, performance or achievements to be materially different from any future results, performance or achievements expressed or implied by the forward-looking statements. The risks and uncertainties that we face are described in our most recent filings with the Securities and Exchange Commission. Except as required by law, we assume no obligation to update these forward-looking statements, even if new information becomes available in the future.

- I. Strategic and Financial Progress – Kevin Hrusovsky – Chairman, CEO
 - i. Q1 Highlights
 - ii. 2019 Goals & Priorities
 - iii. Transforming Medicine: Neurology Momentum Update
- II. Financial Report – Amol Chaubal – CFO
- III. Summary of QTRX Opportunity - Kevin Hrusovsky
- IV. Q&A

Massive Market Opportunity, Disruptive Technology and Continued Growth in Revenues & Publications



Neuro + Oncology



Total QTRX Revenue Growth

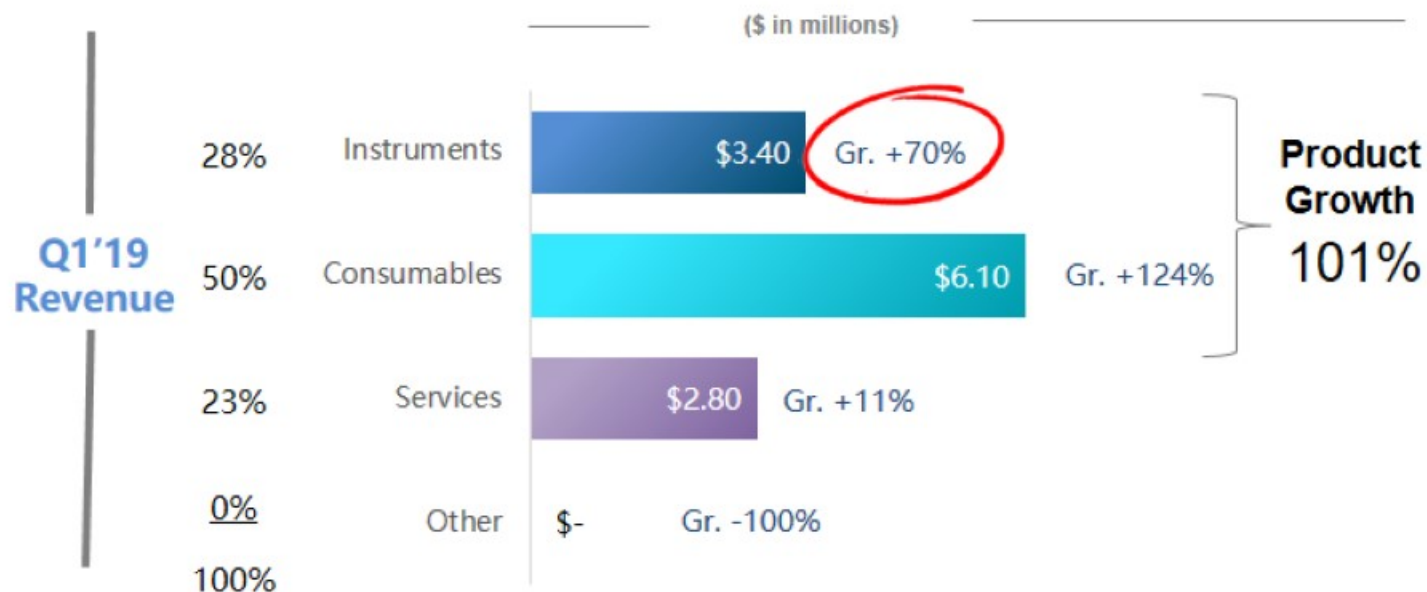


Q1 Highlights

- 64% Revenue growth YOY
- 49% Gross Margin vs 42% PY
- Hired Amol Chaubal, CFO, Mary Ellen Cortizas, Accelerator/CLIA Lab
- Launched SP-X, CorPlex 10 Plex, 9 m-plex assays, 52 onco / immune markers
- 61 new Simoa peer reviewed publications, bringing total > 500
 - sNf-L elevation detected in Alzheimer's 16 years before symptoms
 - LIF biomarker offers promise for blood test for pancreatic cancer

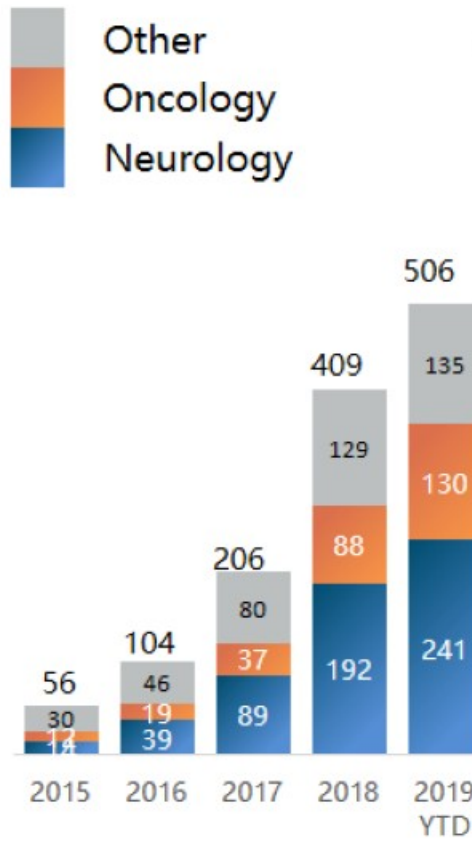
Q1 2019 Growth Led by Consumables and Instruments

- **64% Growth**
- **High margin mix accelerating**
- Consumables growth +124%; GM 48.2%, +450bp
- Total GM 48.7%, +650bp



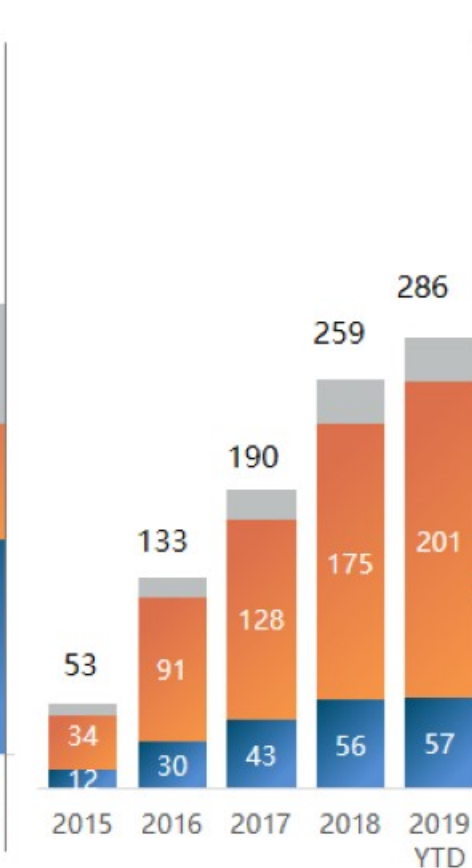
Scientific Research is Driving Brand Awareness, Performance and Utilization

PUBLICATIONS**



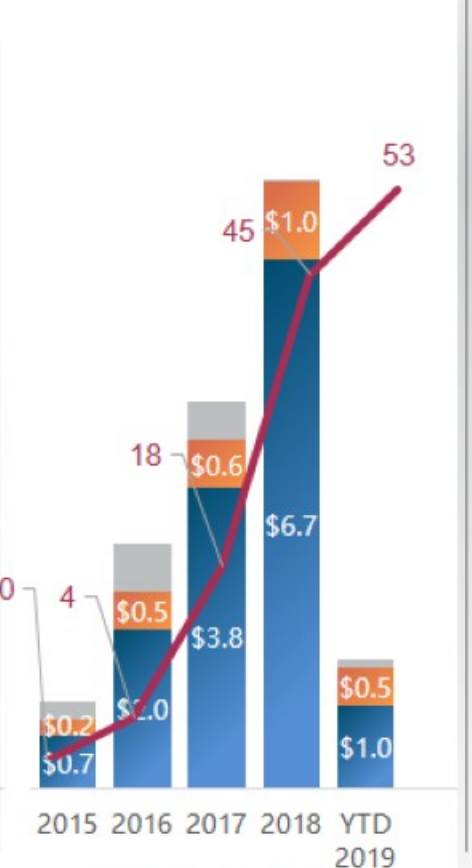
**Cumulative

MARKERS**



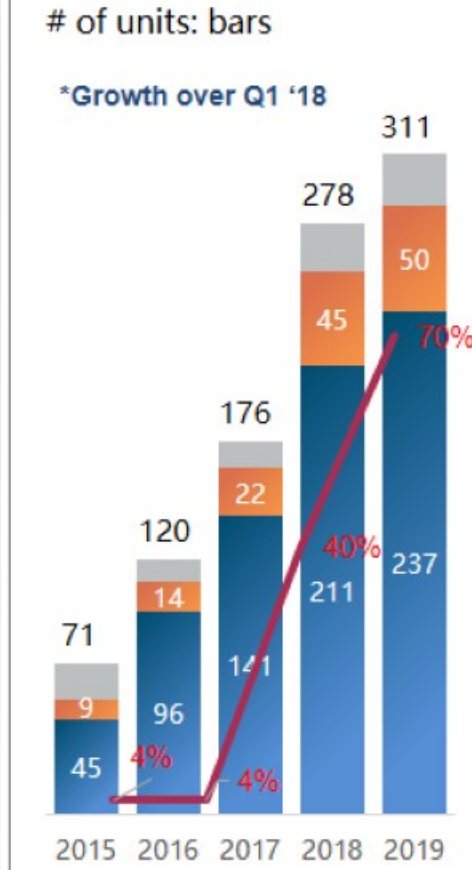
**Cumulative

ACCELERATOR



— Number of drug trial projects

INSTRUMENTS*



— Revenue Growth %

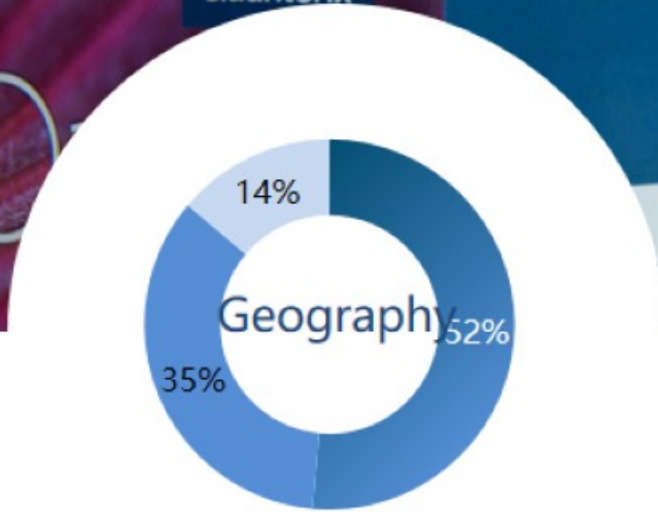
*Instrument segmentation estimated based on consumables sold

CONSUMABLES

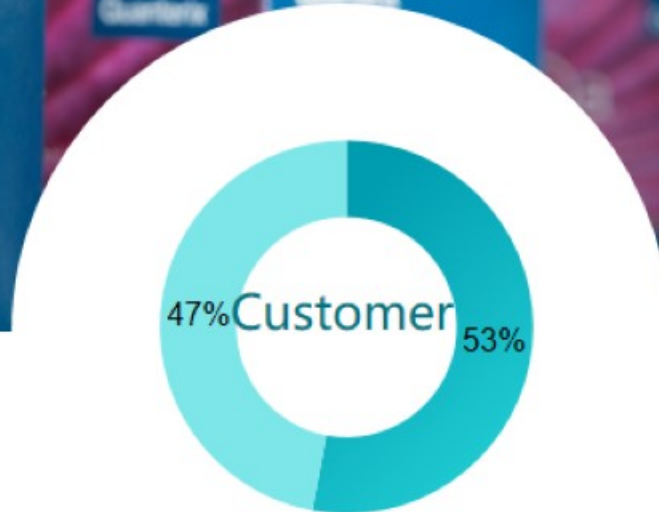


— Consumable growth

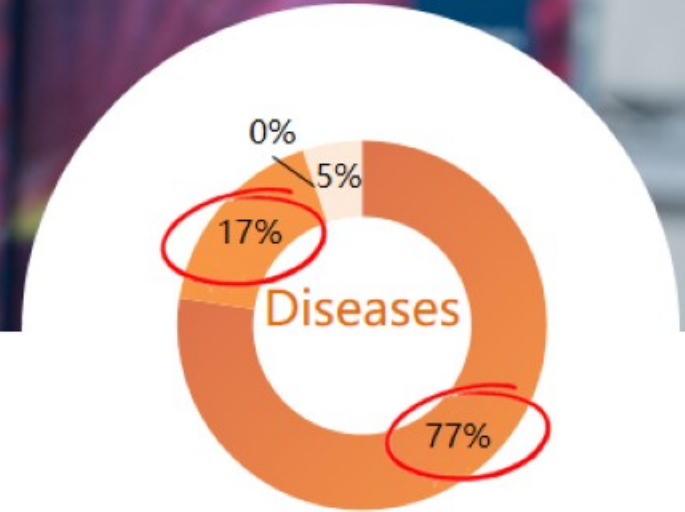
Q1 2019 Growth Stratification



| | <u>Growth</u> |
|--------|---------------|
| NA | +93% |
| Europe | +96% |
| Asia | +161% |



| | <u>Growth</u> |
|--------------------|---------------|
| Pharma/Biotech CRO | +126% |
| Academia | +79% |



| | <u>Growth</u> |
|------------|---------------|
| Neurology | +141% |
| Oncology | +86% |
| Cardiology | -82% |
| ID | +91% |

} 95%

Neurology

High double digit growth

<10% penetrated

Add 25 assays, launch HD-X 4Q'19

Scale globally, LDT/CRO, IT and facilities. Add 50 HC

Oncology

3x Neuro

New entrant

Launch SP-X, CorPlex Assays – 10 plex. 2Q'19

Diagnostics:

50 Drug Trials (neuro/onco)

LDT direct / partner

Explore IVD direct / partner

PPH

Financials:

Growth: 40%

Inst Growth: 25%

Gross Margin: 300 bps

Consumable Utilization 33% of list price or above

Next Generation



- **100x more sensitivity targeted by 2021 YE**
- Lead next frontier of medicine; Protein Translational Modifications

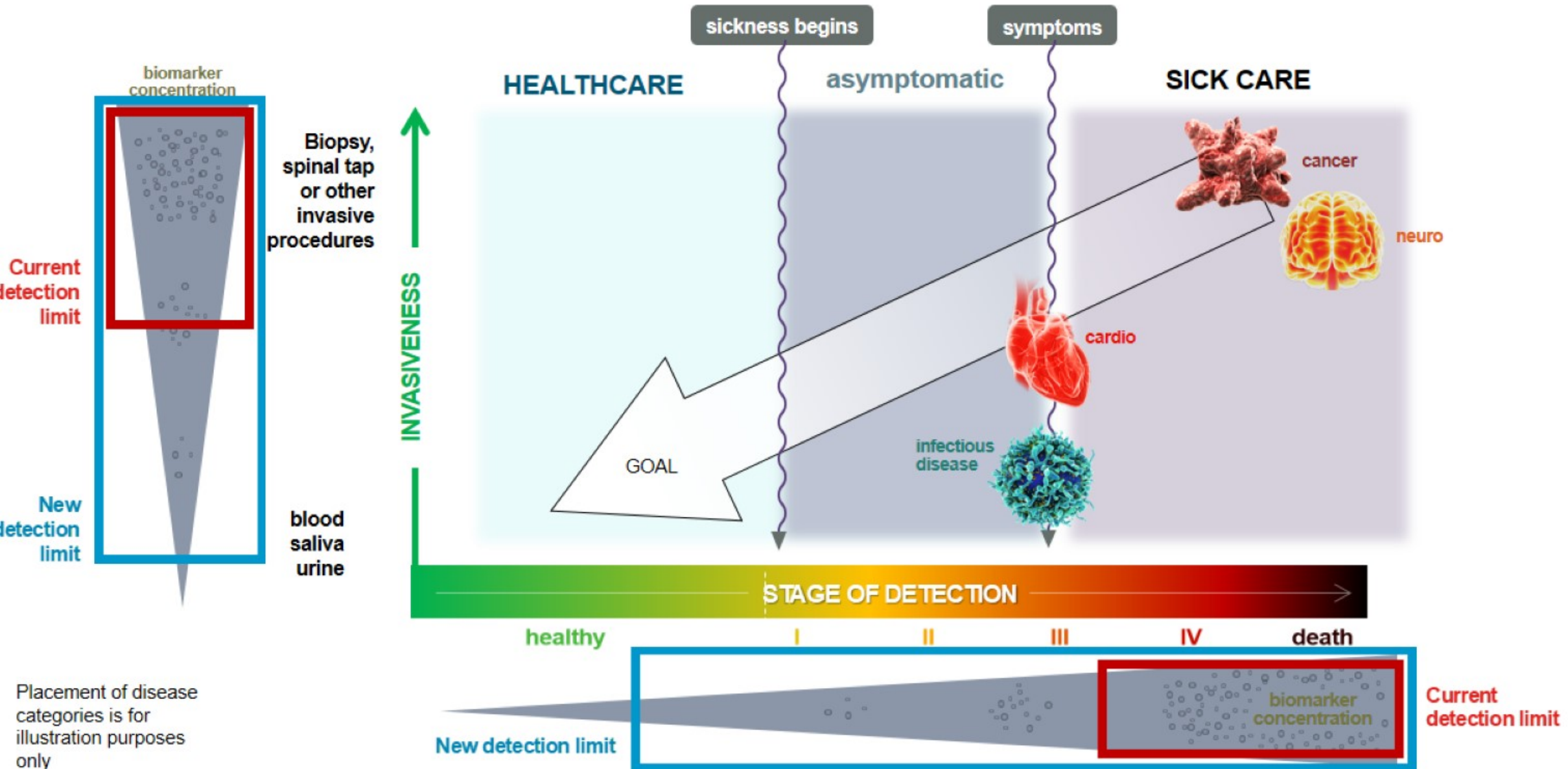
Today: Invasive and Late



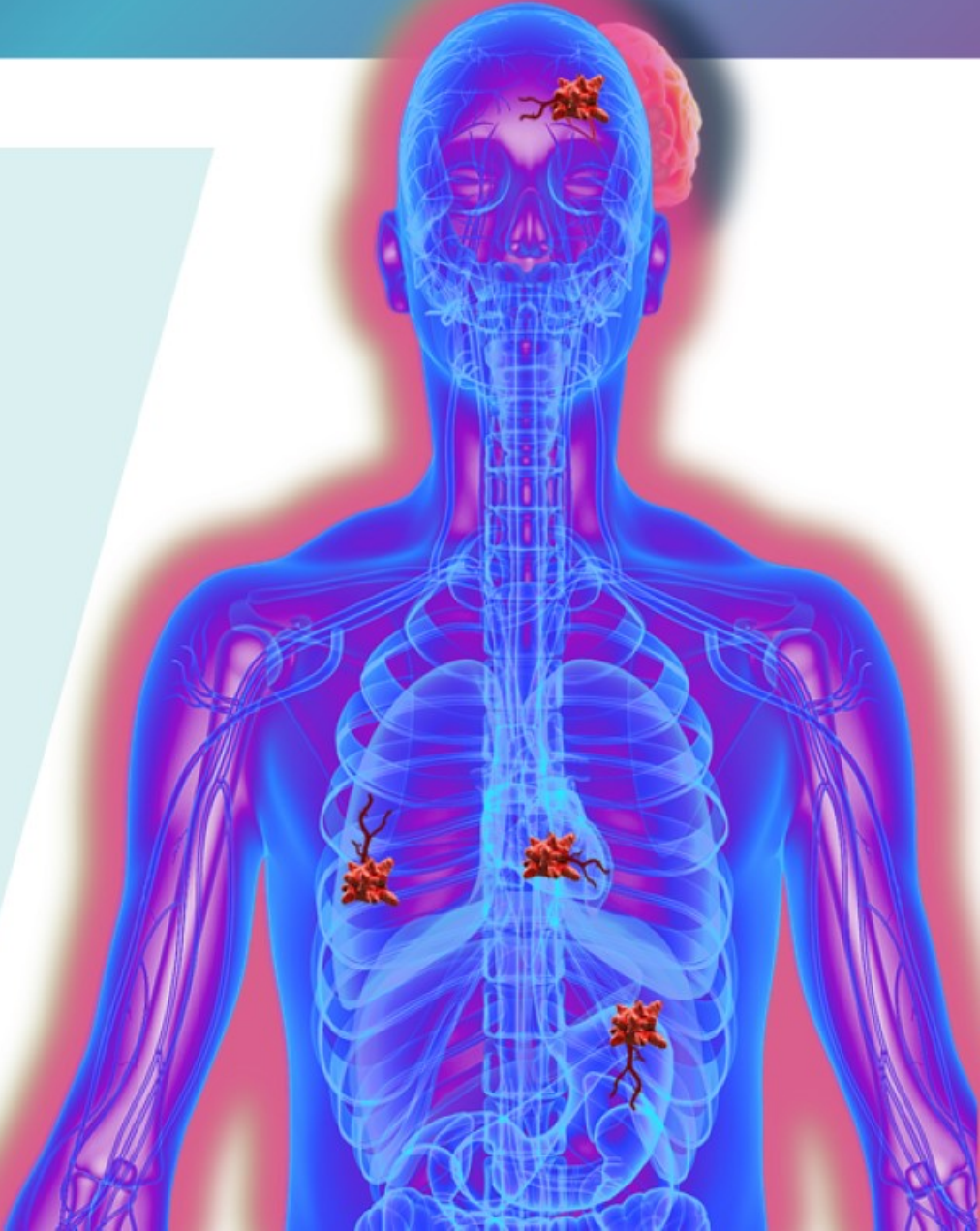
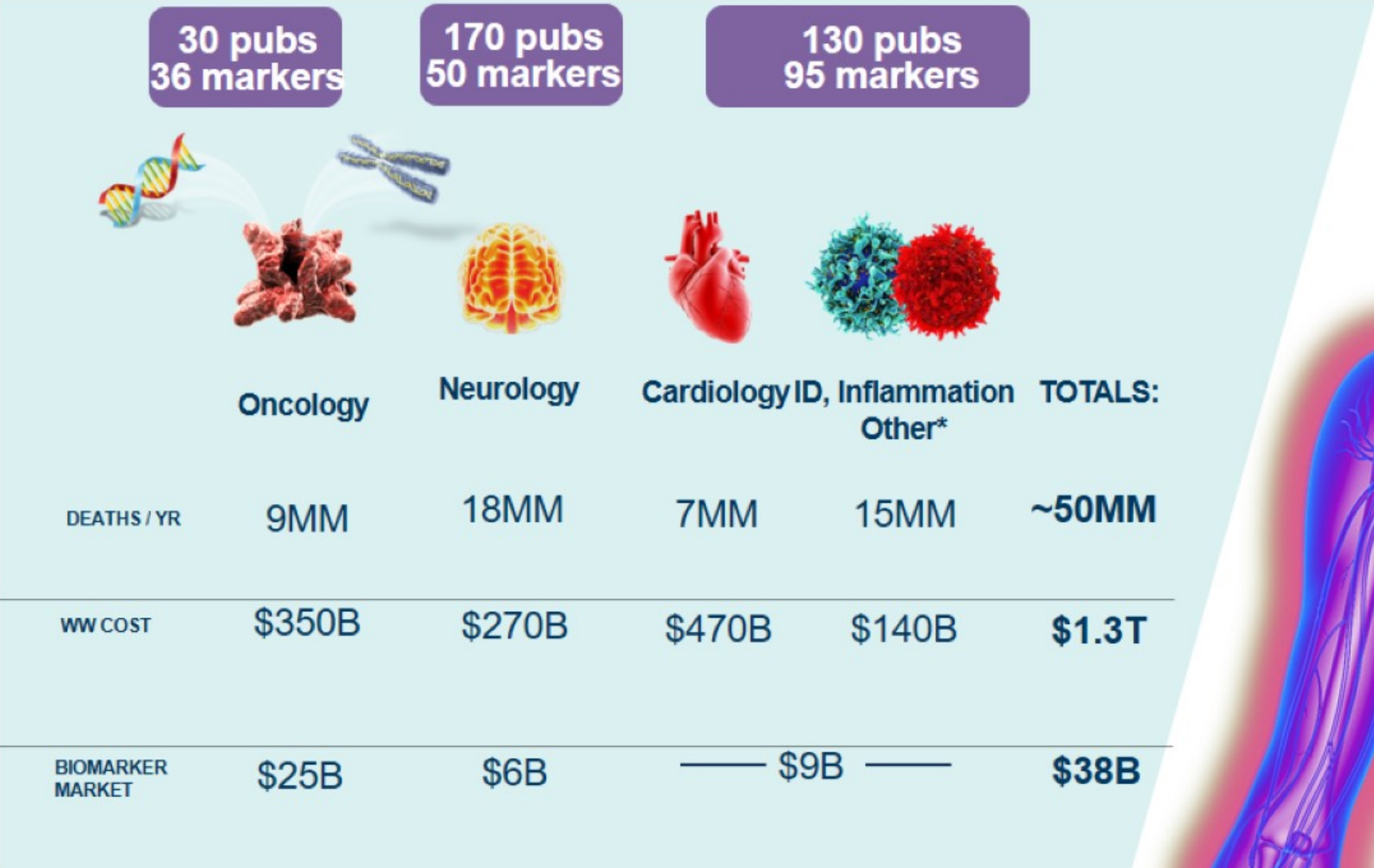
Tomorrow: Non-invasive and Early



Simoa Sees Health to Disease Continuum



BIOMARKER DISRUPTION - Applications



Instruments

Assay kits

Services



HD-1 / HD-X Q4'19

Floor-standing integrated automated system

Assay prep and detection (sample->answer)

400+ publications



SR-X

Benchtop semi-automated assay prep using standardized benchtop devices



SP-X Q2'19

Ultra-sensitive Simoa planar assay technology

Benchtop semi-automated assay prep

Unique multiplex capabilities



Plate based Bead based

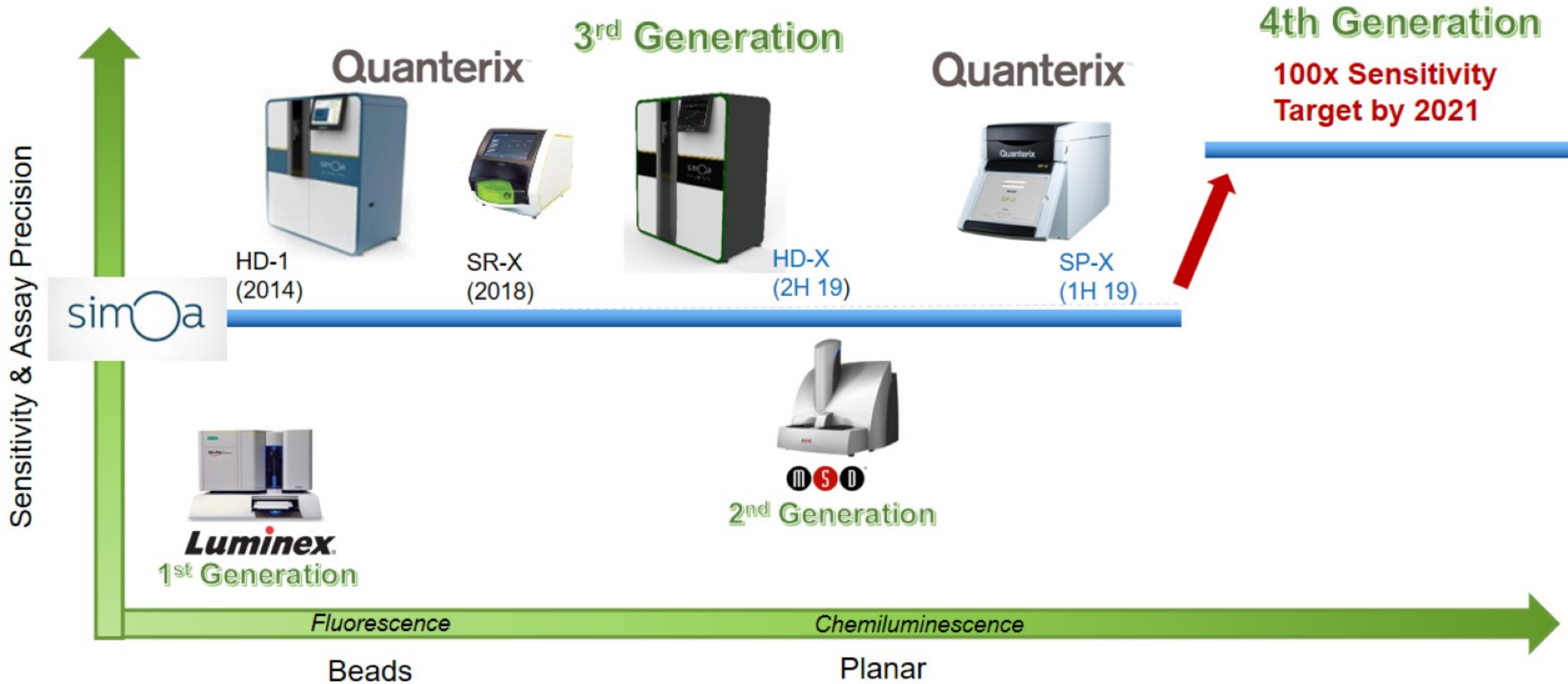
250+ assays for neurology, oncology, cardiology, infectious disease and inflammation

Customization with homebrew kits

Singleplex and multiplex formats



- ✓ Contract research services through Simoa Accelerator Laboratory
- ✓ Sample testing services
- ✓ Custom assay development
- ✓ Custom reagent production and kitting
- ✓ **CLIA and LDT capabilities**



| | Quanterix | | Luminex | MILLIPORE SIGMA Singulex | M S O | protein simple |
|--------------------------|----------------------|----------------------|------------------------|-----------------------------|---------------------------|----------------|
| Technology | Bead-Based Simoa | Planar SP-X Simoa | Bead-based Immunoassay | Erenna Immunoassay | Electrochemi-luminescence | ELLA |
| Sensitivity | Green | Yellow (highlighted) | Red | Green | Red | Red |
| Dynamic Range | Green | Green | Green | Yellow | Green | Yellow |
| Automation & Ease of Use | Green | Yellow (highlighted) | Yellow | Red | Yellow | Green |
| Precision | Green | Green | Red | Red | Yellow | Green |
| Multiplexing | Yellow (highlighted) | Green | Green | Red | Green | Yellow |
| Menu | Yellow (highlighted) | Yellow (highlighted) | Green | Red | Green | Yellow |
| Cost | Yellow (highlighted) | Green | Yellow | Red | Yellow | Yellow |

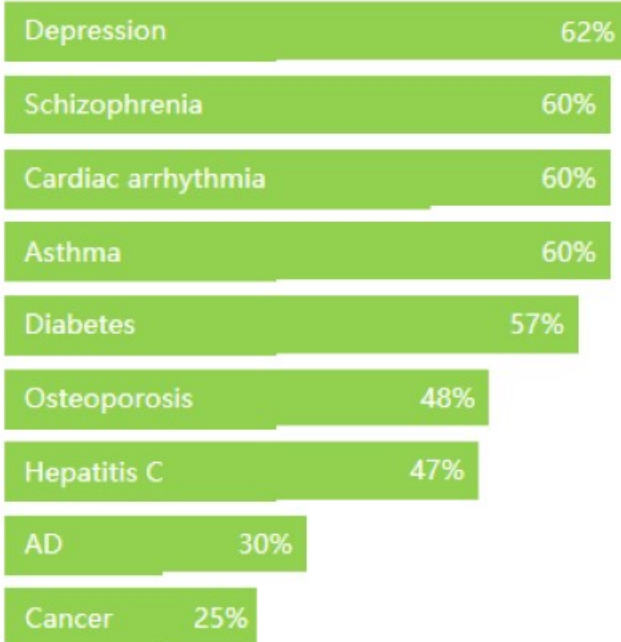
FDA Announces Office of Drug Evaluation Science - ODES

DRUG PERFORMANCE

TOXICITY

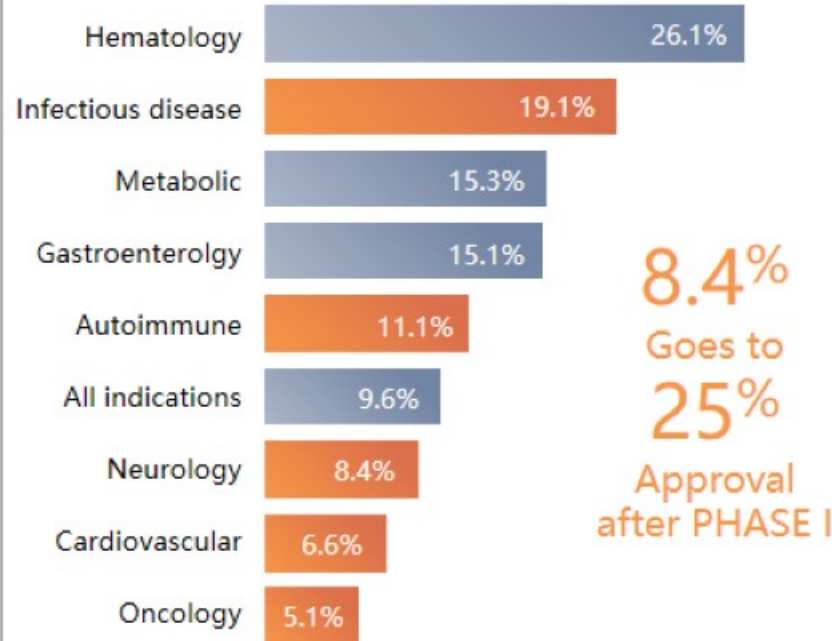
Adverse drug events are a substantial cause of Death in USA

EFFICACY



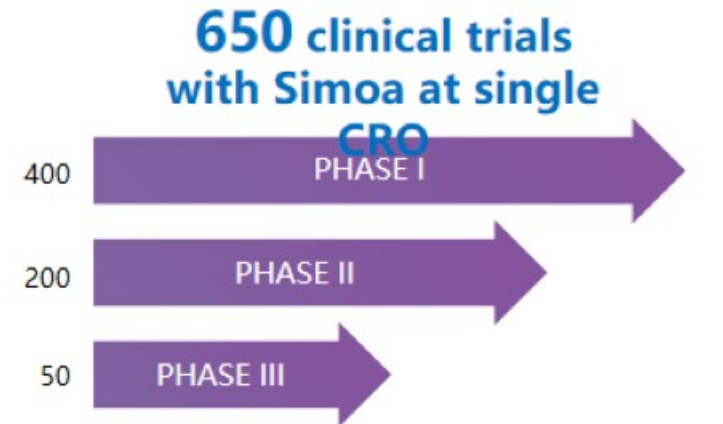
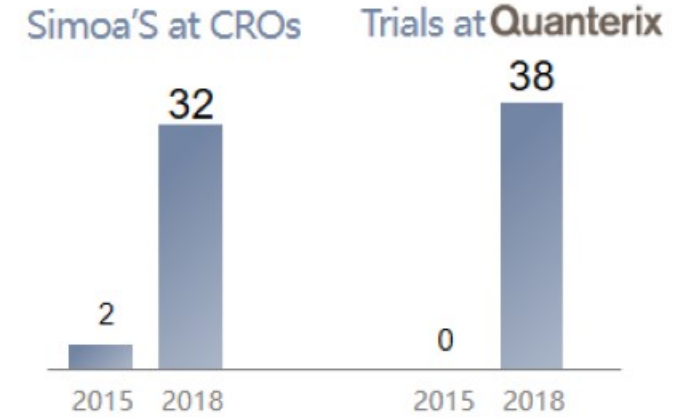
PROBABILITY OF DRUG APPROVAL

300%
increase if biomarkers are used



Probability of phase III approval after Phase 1 approval

VALIDATION OF SIMOA IMPACT



MYRIAD RBM.

Early Alzheimer's Disease: Developing Drugs for Treatment Guidance for Industry

*Contains Nonbinding Recommendations
Draft - Not for Implementation*

1 **Breakthrough Devices Program**
2 **Draft Guidance for Industry and**
3 **Food and Drug Administration Staff**

FDA INITIATIVE...

WHAT IT MEANS FOR QTRX...

Recognizes the criticality of using biomarkers.

Encourages biopharma to identify/incorporate promising biomarkers in trial design



GUIDANCE DOCUMENT
Alzheimer's Disease: Developing Drugs for Treatment Guidance for Industry FEBRUARY 2018
ENDPOINTS FOR STAGE 3,2,1 TRIALS
CDER: BILLY DUNN

FEDERAL REGISTER
The Daily Journal of the United States Government

Program for Parallel Review of Medical Devices

A Notice by the Centers for Medicare & Medicaid Services and the Food and Drug Administration on 10/24/2016

DEPARTMENT OF HEALTH AND HUMAN SERVICES
Centers for Medicare & Medicaid Services
[CMS-3180-N4]
Food and Drug Administration
[Docket No. FDA-2016-N-0308]
Program for Parallel Review of Medical Devices

AGENCY: Food and Drug Administration; Centers for Medicare & Medicaid Services, HHS.
ACTION: Notice.

SUMMARY: The Food and Drug Administration (FDA) and the Centers for Medicare & Medicaid Services (CMS) (the Agencies) are informing the public that the Parallel Review of medical devices pilot program will be fully implemented and extended indefinitely. The Agencies are soliciting nominations from manufacturers of innovative medical devices to participate in the "Program for Parallel Review of Medical Devices." The Parallel Review program is a collaborative effort that is intended to reduce the time between FDA marketing approval or FDA's granting of a de novo request and Medicare coverage decisions through CMS's National Coverage Determination (NCD).

Expedited development, assessment, and review

Expedited review of IVD tests; more probable Class II path. Example: Banyan TBI test (6 mo to clearance).

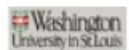
FDA teaming up with CMS to provide parallel review of medical device approval and Medicare coverage.

Potential to secure reimbursement determination simultaneous with FDA clearance

Research Institutions

Biopharma

Other



Vejle Hospital
- part of Lillebælt Hospital



MYRIAD RBM



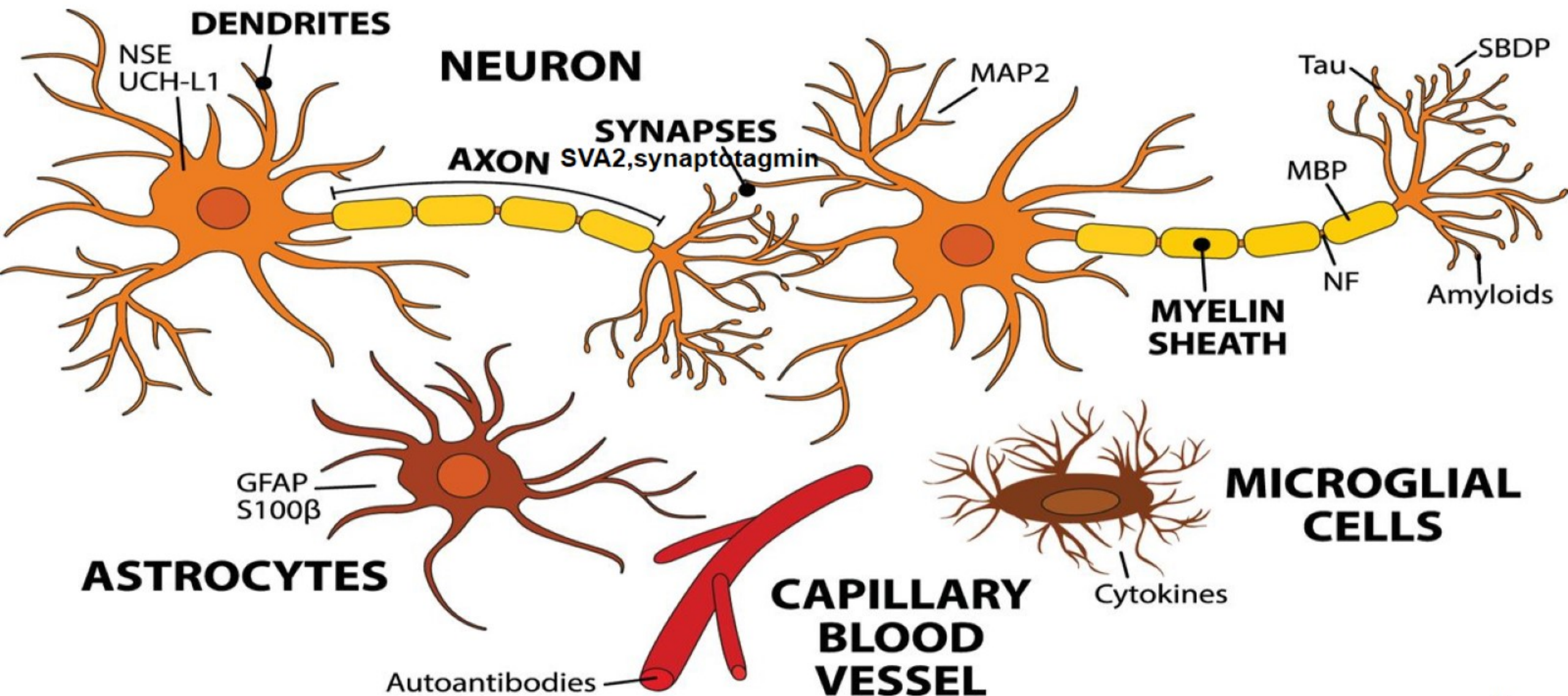
Blue Chip Customers

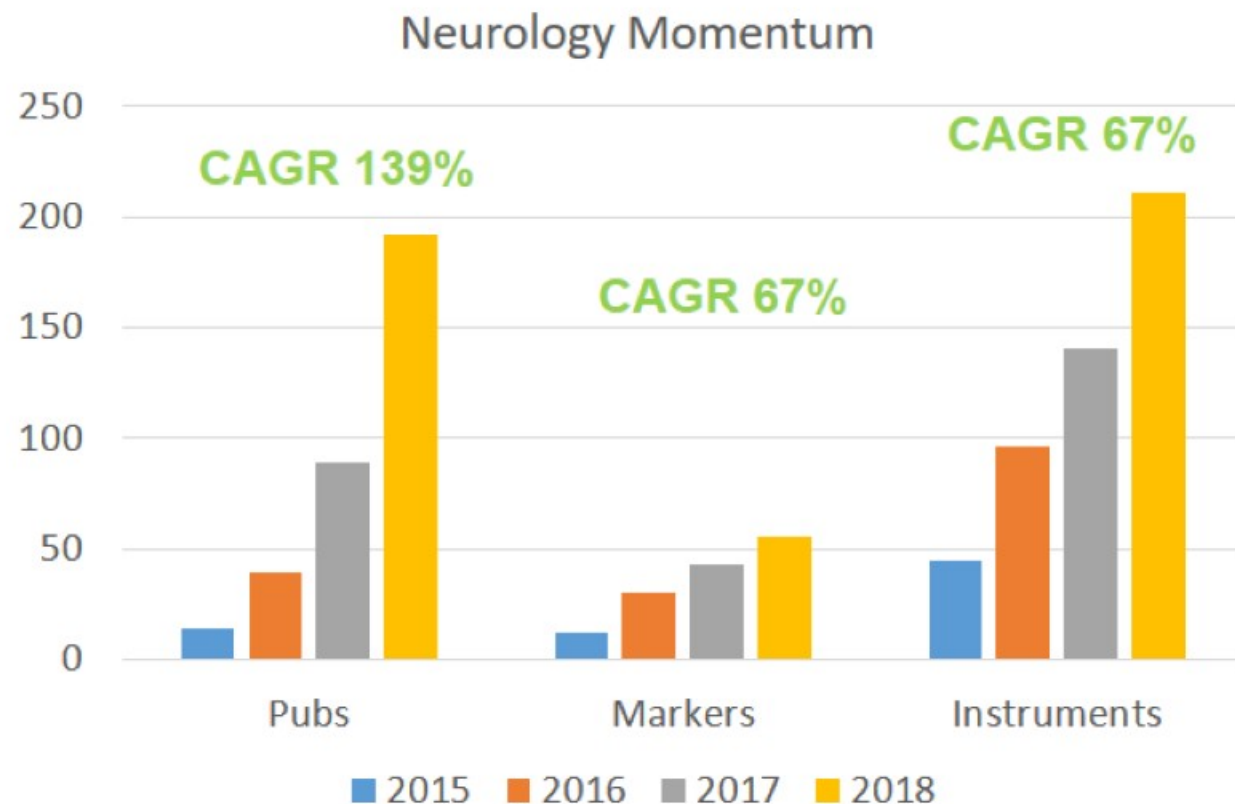
Research Institutions

Biopharma

Other



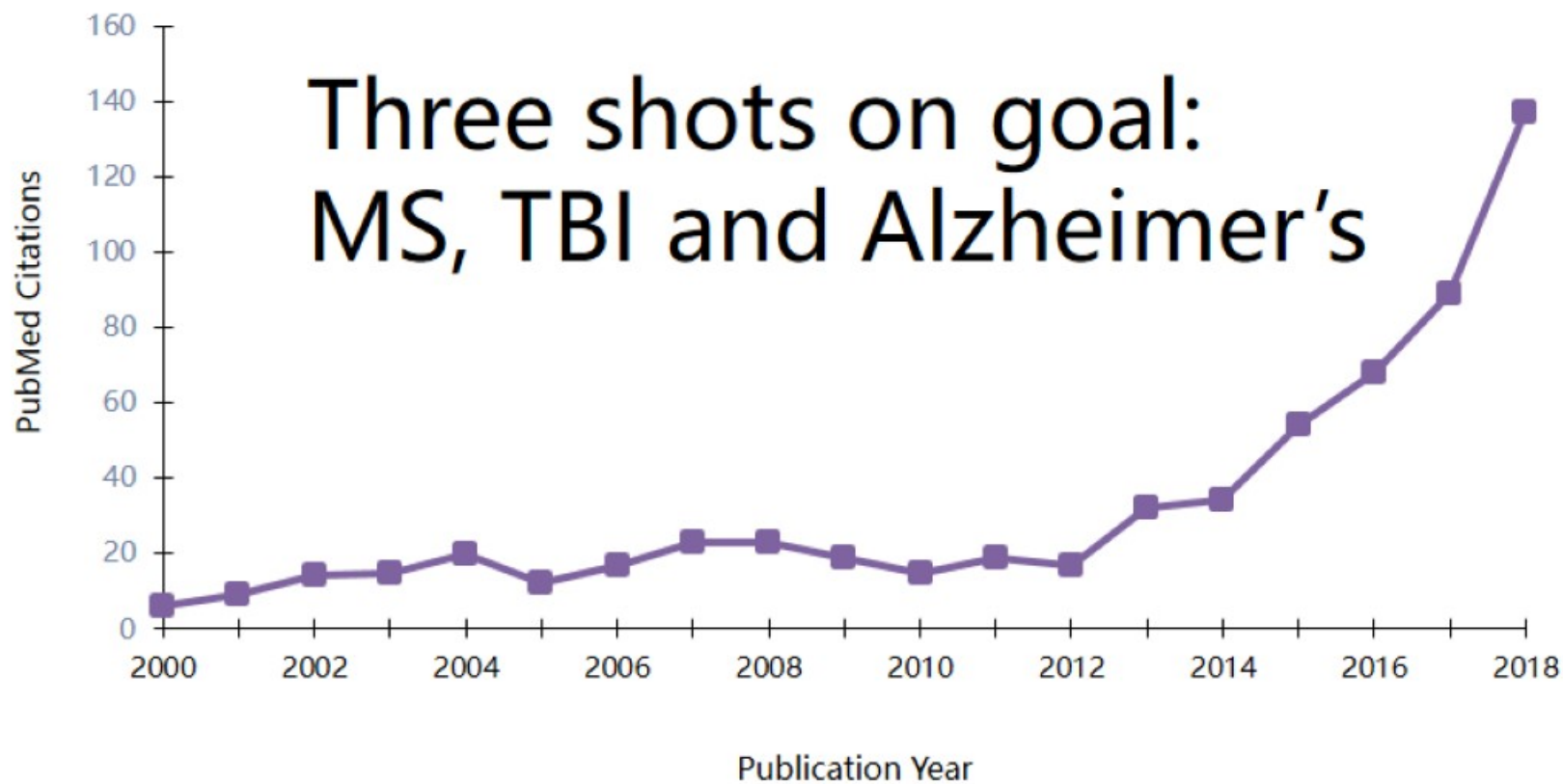




- Strongest **growth in neurology** RUO franchise with **accelerating publications** 2018
- **Nf-L, amyloid beta, tau**, and other markers showing tremendous **potential for key clinical applications** in top peer-reviewed pubs
- Strong network of **world-leading KOLs**
- Inbound **FDA interest** (Nf-L); recent **breakthrough device exemption** for AB40/42 mass spec test for Alzheimer's
- **Media, patient advocacy, and public attention** to promise of blood-based neurology tests (CNN, Forbes, Bloomberg, Washington Post, GMA, etc.)
- **Deep adoption** by leading academic and pharma



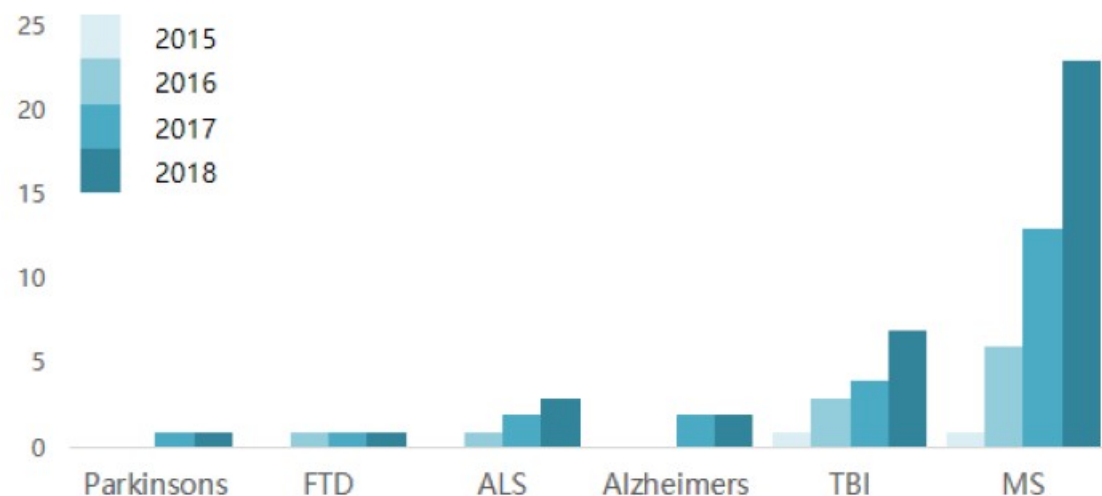
RESEARCH PUBLICATIONS ON NFL



46 active clinical trials using sNf-L



NFL PUBLICATIONS

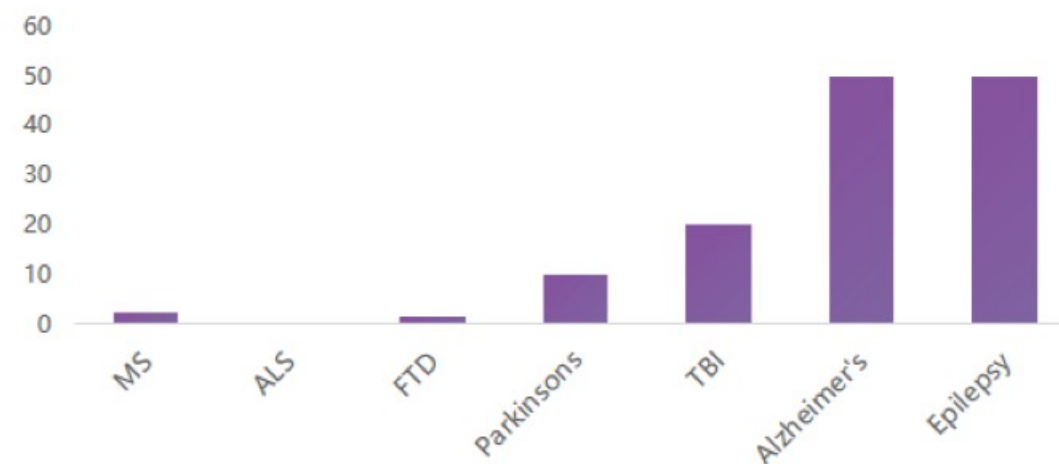


STUDIES CONFIRM NFL CLINICAL UTILITY:

- Disease activity monitoring
- Drug efficacy monitoring
- Relapse/severity prognostic

Majority of published data obtained with Simoa NfL

WW DISEASE INCIDENCE (MILLIONS)



MULTIPLE SCLEROSIS:

- Avg. age of onset: 34 yrs; avg. life expectancy after onset: 30 yrs
- Standard of care: MRI 1-2X/yr
- NfL as MRI replacement: 3.5M tests/yr

Clinical Validation of NfL for MS is a Key Beachhead

Neurology®

JAMA Neurology

Blood neurofilament light chain as a biomarker of MS disease activity and treatment response

Jens Kuhle, MD,* Harald Kropshofer, PhD,* Dieter A. Haering, PhD, Uma Kundu, MPharm, Rolf Meinert, PhD, Christian Barro, MD, Frank Dahlke, MD, Davorka Tomic, PhD, David Leppert, MD, and Ludwig Kappos, MD
Published February 8, 2019

Conclusions:

Blood NfL levels are associated with clinical and MRI-related measures of disease activity and neuroaxonal damage and have prognostic value. Our results support the utility of blood NfL as an easily accessible biomarker of disease evolution and treatment response.

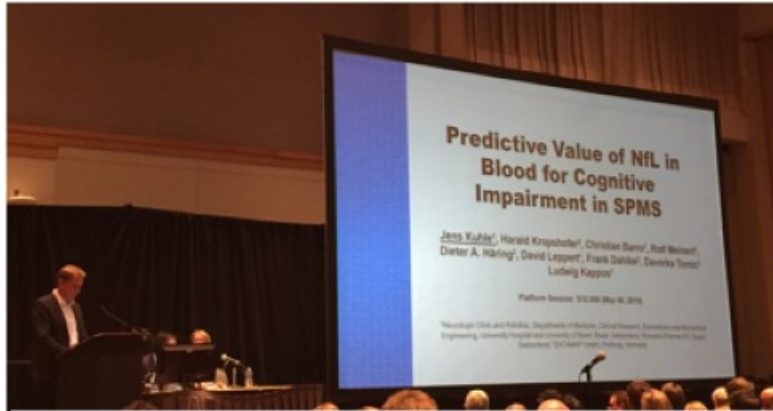
Association Between Longitudinal Plasma Neurofilament Light and Neurodegeneration in Patients With Alzheimer's Disease

Niklas Mattsson, MD, PhD; Nicholas C. Cullen, BSc; Ulf Andreasson, PhD; Henrik Zetterberg, MD, PhD; Kaj Blennow, MD, PhD
JAMA Neurol. doi:10.1001/jamaneurol.2019.0765

Conclusions:

The findings suggest that plasma NfL can be used as a noninvasive biomarker associated with neurodegeneration in patients with AD and may be useful to monitor effects in trials of disease-modifying drugs.

AAN 2019: Explosive momentum of Simoa sNf-L



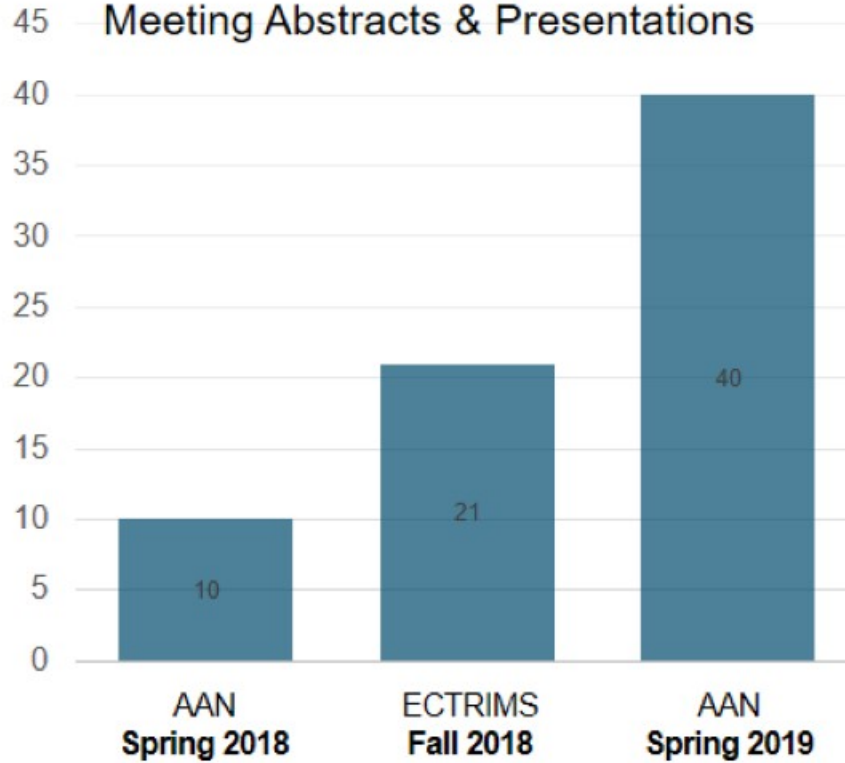
Quanterix' Simoa Technology Powers More Than 85 Percent of Neurofilament Light Biomarker Research to be Unveiled at American Academy of Neurology Annual Meeting

Leading global neurology conference will feature 36 new Simoa-powered studies validating the use of Neurofilament light chain (NFL) as a potential diagnostic and prognostic biomarker for neurodegeneration

May 06, 2019 11:30 AM Eastern Daylight Time



Simoa NFL Meeting Abstracts & Presentations



5 Phase III trials across >1600 MS patients demonstrate clinical utility of Simoa NFL for relapsing remitting MS disease monitoring

Serum Neurofilament Light (NFL) for Disease Prognosis and Treatment Monitoring in Multiple Sclerosis Patients: Toward Implementation Into Clinical Care.

Peter A. Calabresi,¹ Jens Kuhle,² Douglas L. Arnold,³ R. Philip Kinser,⁴ Ludwig Kappos,⁵ Carol M. Singh,⁶ Dipen Sangundekar,⁷ Carl De Moor,⁸ Seth Engle,⁹ Ray Su,¹⁰ Aaron Deykin,¹¹ Elizabeth Fisher,¹² Alfred Sandrock,¹³ Bernd C. Kieseier,¹⁴ Richard A. Rudick,¹⁵ Tatiana Pavesi¹⁶

¹Department of Neurology, Johns Hopkins University School of Medicine, Baltimore, MD, USA; ²Neurologic Clinic and Polyclinic, Departments of Medicine, Biomedicine and Clinical Research, University Hospital Basel, Basel, Switzerland; ³Montreal Neurological Institute, McGill University, Montreal, QC, Canada; ⁴Department of Neurosciences, University of California, San Diego, CA, USA; ⁵Biogen Inc., Cambridge, MA, USA



Natalizumab Reduces Serum Concentrations of Neurofilament Light Chain in Secondary Progressive Multiple Sclerosis Patients From the Phase 3 ASCEND Study

Raju Kapoor,¹ Finn Sellebjerg,² Hans-Peter Hartung,³ Douglas Arnold,⁴ Mark S. Freedman,⁵ Douglas Jeffery,⁶ Aaron Miller,⁷ Keith R. Edwards,⁸ Carol M. Singh,⁹ In Chang,¹⁰ Zhang Ren,¹¹ Dipen Sangundekar,¹² Bing Zhu,¹³ Devangi Kheria,¹⁴ Pei-Ran Ho,¹⁵ Felicia Campbell,¹⁶ Michael Edwards,¹⁷ Elizabeth Fisher,¹⁸ Bernd C. Kieseier,¹⁹ Richard A. Rudick,²⁰ Tatiana Pavesi²¹

Long-term Effect of Fingolimod in Reducing Blood Neurofilament Light Levels in Patients with Relapsing-remitting Multiple Sclerosis

Jeffrey Cohen,¹ Ludwig Kappos,² Nadia Tenenbaum,³ Jackie Han,⁴ Harald Kropshofer,⁵ Devorka Tomić,⁶ Jens Kuhle⁷

¹Cleveland Clinic, ²Neurologic Clinic and Polyclinic, Department of Medicine, Clinical Research, Biomedicine and Biomedical Engineering, University Hospital and University of Basel, ³Novartis Pharmaceuticals Corporation, ⁴Novartis Pharma AG

Objective: To assess the effect of long-term treatment with fingolimod on blood neurofilament light chain (NFL) levels in patients with relapsing-remitting multiple sclerosis (RRMS).

Background: NFL, a cytoskeleton protein, is elevated in blood upon neuroaxonal damage. Blood NFL is a promising biomarker for monitoring disease activity, treatment response, and prognosis in MS.

Design/Methods: This post hoc analysis was based on data from patients who received fingolimod 0.5 mg once daily or placebo/interferon beta-1a (IFN) 30 µg once weekly in pivotal studies (24-month FREEDOMS/12-month TRANSFORMS), and then fingolimod in the open-label LONGTERMS extension study for up to 10 years. The analysis included a subset of patients who had blood NFL assessments at baseline, end of core (EoC) in pivotal studies, and end of study (EoS) in LONGTERMS. Patients were categorized into two groups: a continuous group (n=37) who received fingolimod throughout the studies and a switched group (n=42) who transitioned from placebo/IFN group to fingolimod in the LONGTERMS. NFL was measured using Single Molecule Array (SIMOA™) immunoassay. The geometric mean change in NFL levels from baseline to EoS was analyzed using Wilcoxon signed-rank test.

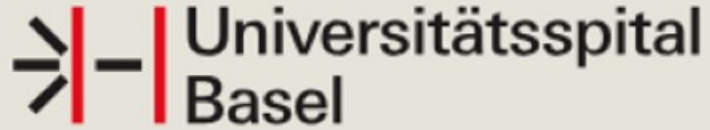
Results: The mean exposure to fingolimod was 3483 days in the continuous group and 2622 days in the switched group. In the continuous group, baseline NFL levels of 23 pg/mL were significantly reduced by approximately 40% at both EoC and EoS (20 pg/mL, P<0.0001 and P=0.0002, respectively). In the switched group, baseline NFL levels of 29 pg/mL were reduced by 15% at EoC (25 pg/mL, P=0.44) and 41% at EoS (17 pg/mL, P<0.0001).

Conclusions: Fingolimod 0.5 mg significantly reduced blood NFL, maintaining its low levels with continuous treatment for up to 10 years. NFL levels were reduced to a lesser extent during treatment with IFN but decreased further with switch to fingolimod, demonstrating the greater impact of highly effective therapy in RRMS.

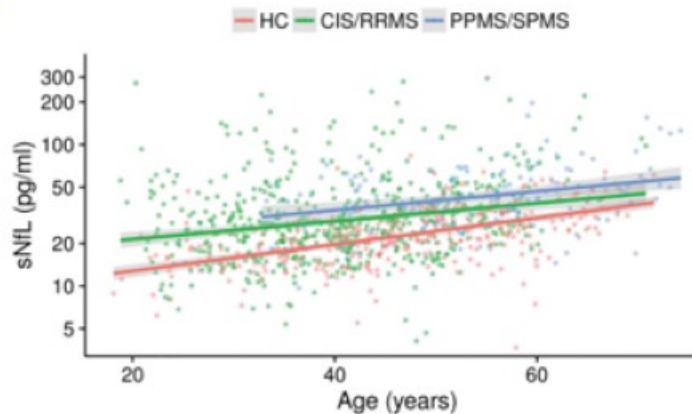
Simoa NfL Multicenter Analytical Validation - 17 sites across Europe and North America

| | Location | Investigator | Study Site |
|----|--------------|---|---|
| 1 | Amsterdam | Dr. Charlotte | Neurochemistry Laboratory, Amsterdam University |
| 2 | Barcelona | Dr. Manuel | Laboratori de Neuroimmunologia Clínica Centre |
| 3 | Basel | Prof. Jens Kuhle, Dr. Zuzanna | Department Biomedicine, Univ Hospital Basel, Switzerland |
| 4 | Dresden | Dr. Katja Akgün Dr. Ziemssen, Tjalf | Neuroimmunological lab, Center of Clinical Neuroscience, Dresden, Germany |
| 5 | Gothenburg | Prof. Kaj Blennow Prof. Henrik | Clinical Neurochemistry Lab, Mölndal Hospital, Mölndal, Sweden |
| 6 | Göttingen | Prof. Wolfgang Dr. Niels Kruse | Institut für Neuropathologie, Universitätsmedizin Göttingen, Germany |
| 7 | London | Dr. Lucia Bianchi Prof. Gavin | Dept of Neuroscience & Trauma Blizard Institute Queen Mary Univ of London, UK |
| 8 | London | Dr. Amanda Prof. Henrik | The DRI Fluid Biomarker Laboratory at University College London, United Kingdom |
| 9 | Mainz | Prof. Stefan Bittner | Klinik für Neurologie, Universitätsmedizin Mainz. |
| 10 | Milan | Dr. Corni Giancarlo Prof. Roberto Furlan | Clinical Neuroimmunology Unit - Institute of Experimental Neurology, Milan, Italy |
| 11 | Montpellier | Prof. Markus Otto Dr. Patrick Oeckl | Hôpital St Eloi, Montpellier, France |
| 12 | Ulm | Prof. Markus Otto Dr. Patrick Oeckl | University of Ulm, Ulm, Germany |
| 13 | Bethesda | Dr. Raturaj Dr. Bibi Bielekova | National Institutes of Health, Bethesda, MD |
| 14 | Ottawa | Dr. Simon Thebault Dr. Freedman, Mark Dr. Booth, Ronald | MS Clinical Ottawa Hospital University of Ottawa, Ottawa ON Canada |
| 15 | Philadelphia | Dr. Marcus Handy Dr. Amit Bar-Or | Perelman School of Medicine, University of Pennsylvania, Department of Neurology, Philadelphia, |
| 16 | Lexington | Kevin Hrusovsky Dr. David Wilson | Quanterix Corp, Lexington MA |
| 17 | Framingham | Dr. Matthew Dr. Martin Kramer | Sanofi Genzyme, Framingham MA |





- Prof. Jens Kuhle, MD, PhD, and Prof. David Leppert, PhD
- 10 collaborating institutions
- 18,000 samples across demographics (2019)
- 7,000-8,000 MS samples (2020-2021)
- Key objective: define what is 'normal' NFL for different ethnicities and age groups
- Partially supported by Quanterix
- Will solidify Simoa NFL as the go-to test for screening and monitoring brain health



Serum Neurofilament Light: A Biomarker of Neuronal Damage in Multiple Sclerosis

Giulio Disanto, MD, PhD,¹ Christian Barro, MD,² Pascal Benkert, PhD,³
Yvonne Naegelin, MD,² Sabine Schädelin, MSc,³ Antonella Giardiello, MD,¹
Chiara Zecca, MD,¹ Kaj Blennow, PhD,⁴ Henrik Zetterberg, PhD,^{4,5}
David Leppert, MD,² Ludwig Kappos, MD,² Claudio Gobbi, MD,¹
Jens Kuhle, MD, PhD,² and the Swiss Multiple Sclerosis Cohort Study Group

Objective: Neurofilament light chains (NFL) are unique to neuronal cells, are shed to the cerebrospinal fluid (CSF), and are detectable at low concentrations in peripheral blood. Various diseases causing neuronal damage have resulted in elevated CSF concentrations. We explored the value of an ultrasensitive single-molecule array (Simoa) serum NFL (sNFL) assay in multiple sclerosis (MS).

Methods: sNFL levels were measured in healthy controls (HC, n = 254) and two independent MS cohorts: (1) cross-sectional with paired serum and CSF samples (n = 142), and (2) longitudinal with repeated serum sampling (n = 246, median follow-up = 3.1 years, interquartile range [IQR] = 2.0–4.0). We assessed their relation to concurrent clinical, imaging, and treatment parameters and to future clinical outcomes.

Results: sNFL levels were higher in both MS cohorts than in HC (p < 0.001). We found a strong association between CSF NFL and sNFL ($\beta = 0.589$, p < 0.001). Patients with either brain or spinal (43.4pg/ml, IQR = 25.2–65.3) or both brain and spinal gadolinium-enhancing lesions (62.5pg/ml, IQR = 42.7–71.4) had higher sNFL than those without (29.6pg/ml, IQR = 20.9–41.8; $\beta = 1.461$, p = 0.005 and $\beta = 1.902$, p = 0.002, respectively). sNFL was independently associated with Expanded Disability Status Scale (EDSS) assessments ($\beta = 1.105$, p < 0.001) and presence of relapses ($\beta = 1.430$, p < 0.001). sNFL levels were lower under disease-modifying treatment ($\beta = 0.818$, p = 0.003). Patients with sNFL levels above the 80th, 90th, 95th, 97.5th, and 99th HC-based percentiles had higher risk of relapses (97.5th percentile: incidence rate ratio = 1.94, 95% confidence interval [CI] = 1.21–3.10, p = 0.006) and EDSS worsening (97.5th percentile: OR = 2.41, 95% CI = 1.07–5.42, p = 0.034).

Interpretation: These results support the value of sNFL as a sensitive and clinically meaningful blood biomarker to monitor tissue damage and the effects of therapies in MS.

ANN NEUROL 2017;81:857–870

“There is an urgent unmet need for reliable biomarkers of neurodegeneration. NFL addresses this need.”

- Prof Jens Kuhle

Blood Test May Detect Alzheimer's In Patients 16 Years Before Symptoms Appear

CNN Health • Food • Fitness • Wellness • Parenting • Live Longer

Blood test could detect Alzheimer's up to 16 years before symptoms begin, study says

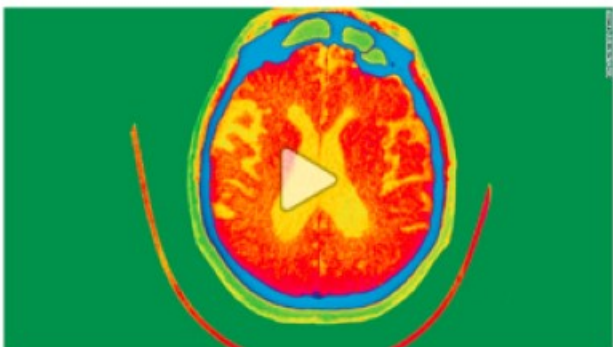
By Nina Aramova, CNN
Updated 2:45 PM ET, Tue January 22, 2019



More from CNN

Passengers spend 14 hours stuck on grounded flight

LFC's Allen Crowder Has 'No Hard Feelings' Toward Greg Hardy...



How to prevent Alzheimer's disease 03:33

(CNN) — A simple blood test could predict if a patient will develop Alzheimer's disease up to 16 years before symptoms begin, a new study finds.

By measuring changes in the levels of a protein in the blood, called neurofilament light chain (NFL), researchers believe any rise in levels of the protein could be an early sign of the disease, according to the study published Monday in the journal Nature Medicine.

NFL is a "marker in the blood which gives an indication of nerve cell loss in the brain," explained lead researcher Mathias Jucker, professor of cell biology of neurological diseases at the German Center for Neurodegenerative Diseases. "The more neurofilament you have in the blood, the more brain damage you have," he said.

nature
medicine

LETTERS

<https://doi.org/10.1038/s41591-018-0304-3>

Serum neurofilament dynamics predicts neurodegeneration and clinical progression in presymptomatic Alzheimer's disease

Oliver Preische^{1,2,3}, Stephanie A. Schultz^{1,2,3}, Anja Apel^{1,2,3}, Jens Kuhle⁴, Stephan A. Kaeser^{1,2}, Christian Barro⁵, Susanne Gräber⁶, Elke Kuder-Buletta⁷, Christian LaFougere⁸, Christoph Laske^{1,3}, Jonathan Vögler^{1,4}, Johannes Levin^{1,4}, Colin L. Masters⁹, Ralph Martins¹⁰, Peter R. Schofield^{11,12}, Martin N. Rossor¹³, Neill R. Graff-Radford¹⁴, Stephen Salloway¹⁵, Bernardino Ghetti¹⁶, John M. Ringman¹⁷, James M. Noble¹⁷, Jasmeer Chhatwal¹⁸, Alison M. Goate¹⁹, Tammie L. S. Benzinger²⁰, John C. Morris²¹, Randall J. Bateman²², Guoqiao Wang²³, Anne M. Fagan²⁴, Eric M. McDade²⁵, Brian A. Gordon²⁶, Mathias Jucker^{1,2,3} and Dominantly Inherited Alzheimer Network²⁰

Neurofilament light chain (NFL) is a promising fluid biomarker of disease progression for various cerebral proteopathies. Here we leverage the unique characteristics of the Dominantly Inherited Alzheimer Network and ultrasensitive immunoassay technology to demonstrate that NFL levels in the cerebrospinal fluid (n = 187) and serum (n = 405) are correlated with one another and are elevated at the presymptomatic stages of familial Alzheimer's disease. Longitudinal, within-person analysis of serum NFL dynamics (n = 196) confirmed this elevation and further revealed that the rate of change of serum NFL could discriminate mutation carriers from non-mutation carriers almost a decade earlier than cross-sectional absolute NFL levels (that is, 16.2 versus 6.8 years before the estimated symptom onset). Serum NFL rate of change peaked in participants converting from the presymptomatic to the symptomatic stage and was associated with cortical thinning assessed by magnetic resonance imaging, but less so with amyloid- β deposition or glucose metabolism (assessed by positron emission tomography). Serum NFL was predictive for both the rate of cortical thinning and cognitive changes assessed by the Mini-Mental State Examination and Logical Memory test.

Thus, NFL dynamics in serum predict disease progression and brain neurodegeneration at the early presymptomatic stages of familial Alzheimer's disease, which supports its potential utility as a clinically useful biomarker.

In most neurodegenerative diseases, brain changes manifest many years before clinical symptoms become apparent. In Alzheimer's disease, presymptomatic changes in the brain include cortical thinning and neuropathological depositions containing amyloid- β and tau. These pathological changes can be assessed by magnetic resonance imaging (MRI), positron emission tomography (PET), and measurement of amyloid- β and tau protein levels in the cerebrospinal fluid (CSF)¹. However, CSF collection is invasive and imaging modalities are expensive; therefore, they are not well suited to routine clinical practice. Blood biomarkers for the presymptomatic phase of Alzheimer's disease are largely lacking, although recent progress in the analysis of amyloid- β , tau, and neurofilament light chain (NFL) in the blood have been reported^{2–5}.

NFL is a component of the axonal cytoskeleton and is primarily expressed in large-caliber myelinated axons^{6,7}. Changes of NFL in bodily fluids have been linked to brain damage and brain atrophy in mouse models and multiple neurological disorders including

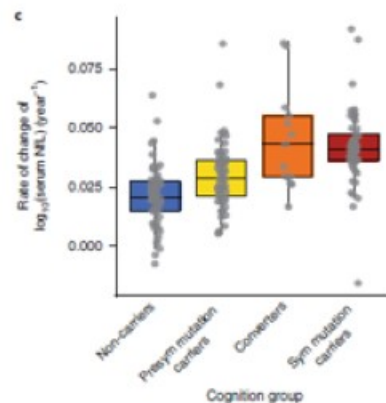


Fig. 3 | Rate of change per year in serum NFL in mutation carriers mirrors rate of change in cortical thinning. a. Relationship between estimated annual rate of change in serum NFL and estimated annual rate of change in precuneus cortical thickness for non-carriers, presymptomatic (Presym) mutation carriers, and symptomatic (Sym) mutation carriers (including converters to the symptomatic phase, see Fig. 2c). Results from LMEMs revealed

¹German Center for Neurodegenerative Diseases (DZNE), Tübingen, Germany. ²Department of Cellular Neurology, Hertie Institute for Clinical Brain Research, and Department of Psychiatry and Psychotherapy, University of Tübingen, Tübingen, Germany. ³Department of Neurology, Department of Radiology, and Division of Biostatistics, Washington University School of Medicine, St. Louis, MO, USA. ⁴Neurologic Clinic and Policlinic, Departments of Medicine, Biomedicine and Clinical Research, University Hospital Basel, University of Basel, Basel, Switzerland. ⁵German Center for Neurodegeneration

University, Providence, RI, USA. ⁶Indiana Alzheimer Disease Center and Department of Pathology and Laboratory Medicine, Indiana University School of Medicine, Indianapolis, IN, USA. ⁷Department of Neurology, Keck School of Medicine at USC, Los Angeles, CA, USA. ⁸Taub Institute for Research on Alzheimer's Disease and the Aging Brain, Department of Neurology, Columbia University Medical Center, New York, NY, USA. ⁹Department of Neurology, Massachusetts General Hospital, Harvard Medical School, Boston, MA, USA. ¹⁰Department of Neuroscience, Icahn School of Medicine at Mount Sinai, New York, NY, USA. ¹¹A full list of members and affiliations appears at the end of the paper. ¹²These authors contributed equally: Oliver Preische, Stephanie A. Schultz, Anja Apel. *e-mail: mathias.jucker@uni-tuebingen.de

Simoa: NfL predicts clinical progression in pre-symptomatic Alzheimer's disease 16 years before onset

Quanterix
The Science of Precision Health

**nature
medicine**

Oliver Preische^{1,2,21}, Stephanie A. Schultz^{3,21}, Anja Apel^{1,2,21}, Jens Kuhle⁴, Stephan A. Kaeser^{1,2}, Christian Barro⁴, Susanne Gräber¹, Elke Kuder-Buletta¹, Christian LaFougere¹, Christoph Laske^{1,2}, Jonathan Vögler^{5,6}, Johannes Levin^{5,6}, Colin L. Masters⁷, Ralph Martins^{8,9}, Peter R. Schofield^{10,11}, Martin N. Rossor¹², Neill R. Graff-Radford¹³, Stephen Salloway¹⁴, Bernardino Ghetti¹⁵, John M. Ringman¹⁶, James M. Noble¹⁷, Jasmeer Chhatwal¹⁸, Alison M. Goate¹⁹, Tammie L. S. Benzinger³, John C. Morris³, Randall J. Bateman³, Guoqiao Wang³, Anne M. Fagan³, Eric M. McDade³, Brian A. Gordon³, Mathias Jucker^{1,2*} and Dominantly Inherited Alzheimer Network²⁰



BREAKING NEWS

NEW BLOOD TEST CAN PREDICT ALZHEIMER'S DISEASE

**LIVE
CNN**

...ED ONLY BY HIS DESIRE TO BRING A SUPER BOWL CHAMPION

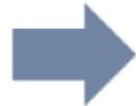
2:12 PM ET

Digital Biomarkers Disruption Paradigm: Alzheimer's Disease Opportunity

Today



Alzheimer's disease not diagnosed until symptoms



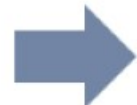
Imaging expensive and often not covered



Therapies for later stage disease have limited effectiveness

Tomorrow

Detect or screen in annual blood test



Follow on Image or blood test to diagnose



Therapy delivered sooner with less dosing / toxicity. Blood test monitors progression

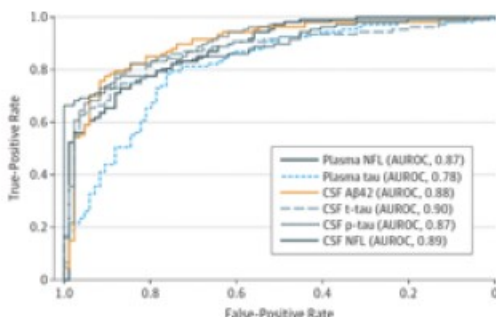


Aducanumab from Biogen

Taking Aim at Alzheimer's Disease with Simple Blood Test

JAMA Neurology | Original Investigation
Association of Plasma Neurofilament Light With Neurodegeneration in Patients With Alzheimer Disease

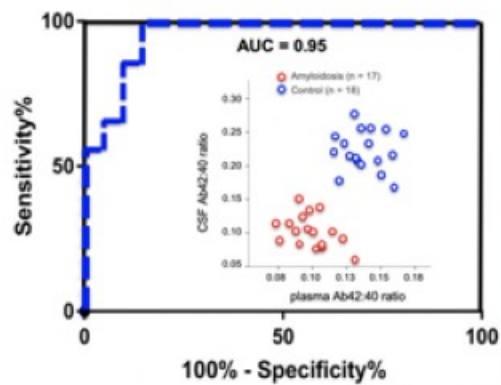
2018



Plasma NfL exhibits high diagnostic accuracy for Alzheimer's Disease (AUC 0.87)

With Sudden Progress, Blood Aβ Rivals PET at Detecting Amyloid

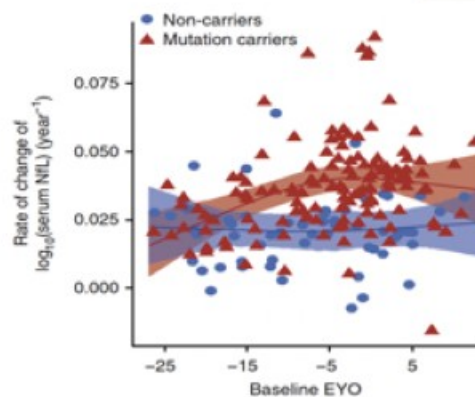
2018



Plasma Aβ42/Aβ40 ratio exhibits high diagnostic accuracy for amyloid positive patients (AUC 0.95)

Serum neurofilament dynamics predicts neurodegeneration and clinical progression in presymptomatic Alzheimer's disease

2019



Serum NfL predicts clinical progression in pre-symptomatic Alzheimer's disease 16 years before onset.

NEW BLOOD TEST CAN PREDICT ALZHEIMER'S DISEASE



Every **65 SECONDS** SOMEONE IN THE UNITED STATES DEVELOPS THE DISEASE
Every **3 SECONDS** SOMEONE IN THE WORLD DEVELOPS THE DISEASE

5.8 MILLION AMERICANS ARE LIVING WITH ALZHEIMER'S. BY 2050, THIS NUMBER IS PROJECTED TO RISE TO NEARLY 14 MILLION.

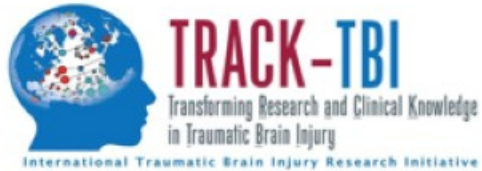
44 MILLION PEOPLE ARE LIVING WITH ALZHEIMER'S WORLDWIDE

LETTERS
https://doi.org/10.1001/ama.2018.0283

Serum neurofilament dynamics predicts neurodegeneration and clinical progression in presymptomatic Alzheimer's disease

Oliver Preische^{1,2,3}, Stephanie A. Schultz^{1,2,3}, Anja Apel^{1,2,3}, Jens Kuhle^{1,2,3}, Stephan A. Kaeser^{1,2,3}, Christian Barro^{1,2,3}, Susanna Graber^{1,2,3}, Elke Kuder-Buletta^{1,2,3}, Christian Lafougere^{1,2,3}, Christoph Laska^{1,2,3}, Jonathan Vogtlein^{1,2,3}, Johannes Levin^{1,2,3}, Colin L. Masters^{1,2,3}, Ralph Martins^{1,2,3}, Peter R. Schofield^{1,2,3}, Martin N. Rossor^{1,2,3}, Neil R. Graff-Radford^{1,2,3}, Stephan Salloway^{1,2,3}, Bernardino Ghetti^{1,2,3}, John M. Ringman^{1,2,3}, James M. Noble^{1,2,3}, Jasmeer Chhatwal^{1,2,3}, Alison M. Goate^{1,2,3}, Tammie L. S. Benzinger^{1,2,3}, John C. Morris^{1,2,3}, Randall J. Bateman^{1,2,3}, Guoqiao Wang^{1,2,3}, Anne M. Fagan^{1,2,3}, Eric M. McDade^{1,2,3}, Brian A. Gordon^{1,2,3}, Mathias Jucker^{1,2,3} and Dominant Inherited Alzheimer Network^{1,2,3}

Neurofilament light chain (NfL) is a promising fluid biomarker of disease progression for various cerebral pathologies. Here we leverage the unique characteristics of the Dominant Inherited Alzheimer Network and ultrasensitive immunoassay technology to demonstrate that NfL levels in the cerebrospinal fluid (CSF) and serum (s-NfL) are correlated with one another and are elevated at the presymptomatic stages of familial Alzheimer's disease. Longitudinal, within-person analysis of serum NfL dynamics (n = 198) confirmed this elevation and further revealed that the rate of change of serum NfL could discriminate mutation carriers from non-mutation carriers almost a decade earlier than cross-sectional amyloid β (Aβ) levels (that is, 16.2 versus 6.8 years before the estimated symptom onset). Serum NfL rate of change peaked in participants converting from the presymptomatic to the symptomatic stage and was associated with cortical thinning assessed by magnetic resonance imaging, but less so with amyloid β deposition or glucose metabolism (assessed by positron emission tomography). Serum NfL was predictive for both the rate of cortical thinning and cognitive changes assessed by the Mini-Mental State Examination and Logical Memory test. Thus, NfL dynamics in serum predict disease progression and brain neurodegeneration at the early presymptomatic stages of familial Alzheimer's disease, which supports its potential utility as a clinically useful biomarker.



Geoff Manley MD PhD
UCSF, TRACK-TBI



Prospective TBI trial comparing Simoa NFL to imaging and outcome measures with generation of large normative database

- Dr. Ramon Diaz-Arristia, principle investigator; TRACK-TBI P.I. Dr. Geoff Manley
- TRACK-TBI: 18 collaborating US sites, 3,000 longitudinal brain injury patients
- Canadian Health Measures Survey / Statistics Canada, 6,000 healthy subjects
- Framingham Heart Study OMNI cohort – ethnic diversity, 900 subjects
- Supported in part by Quanterix

Published Ahead of Print on April 13, 2018 as 10.1212/WNL.0000000000005518

ARTICLE IN PRESS

Neurofilament light and tau as blood biomarkers for sports-related concussion

Pashun Shafiq, MD, PhD, Veeravrat Tegner, MD, PhD, Niklas Marklund, MD, PhD, Kaj Blennow, MD, PhD, and Henrik Zetterberg, MD, PhD

Correspondence: Dr. Shafiq, pashun.shafiq@qu.com

Neurology® 2018;01:9. doi:10.1212/WNL.0000000000005518

Abstract

Objective
To compare neurofilament light (NFL) and tau as blood-based biomarkers for acute sports-related concussion (SRC) and determine whether their concentrations at different time points after the injury are associated with prolonged time to return to play (RTP).

Methods
A total of 288 professional hockey players were followed longitudinally from September 1, 2012, to April 30, 2015. Data collection and biomarker analyses were conducted between 2015 and 2017. Associations were tested between blood concentrations of NFL and tau, and RTP time. Serum concentrations of S100B and neuron-specific enolase (NSE) were also measured for comparison.

Results
Of 288 players, 105 sustained an SRC. Of those, 87 underwent blood sampling 1, 12, 36, and 144 hours after SRC and at the RTP time point. Serum NFL concentrations 1, 12, 36, and 144 hours after SRC were related to prolonged RTP time, and could separate players with RTP >10 days from those with RTP ≤10 days (area under the receiver operating characteristic curve [AUROC] 0.82). Also, serum NFL 144 hours after SRC discriminated players who resigned from the game due to persistent postconcussion symptoms (PCS) from those who returned to play (AUROC 0.89). Plasma tau 1 hour after SRC was related to RTP but less strongly than NFL, while S100B and NSE showed no such associations.

Conclusion
Serum NFL outperformed tau, S100B, and NSE as a biomarker for SRC. From a clinical standpoint, serum NFL may be useful to identify individuals at risk of prolonged PCS, and may aid in biomarker-informed decisions with regard to when RTP should be considered.

Ultrasensitive LIF Assay Revealed as Promising Blood Biomarker Drug Target for Pancreatic Cancer

Science News

from research organizations

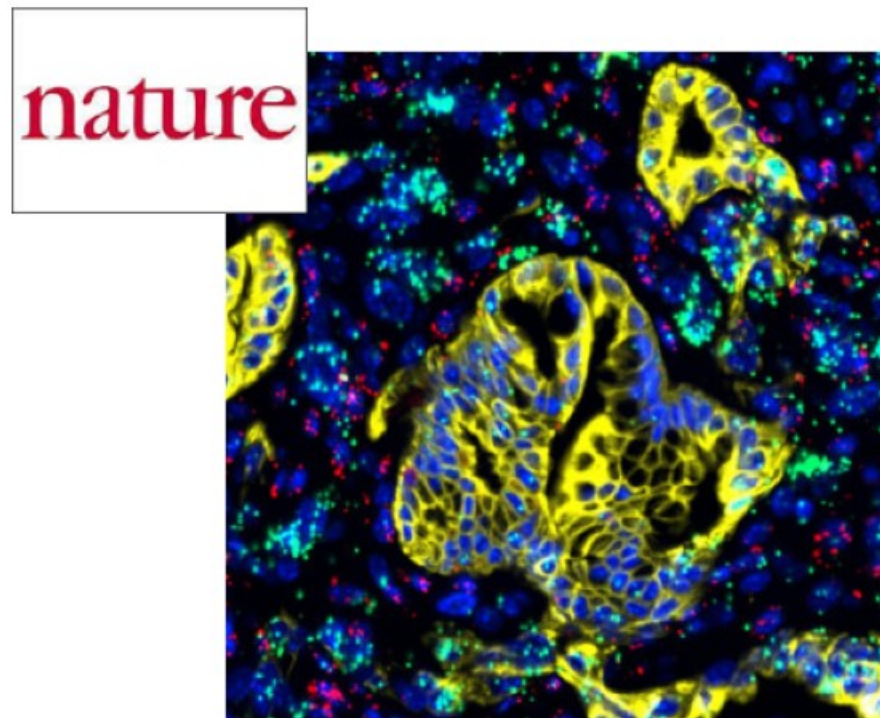
New study targets Achilles' heel of pancreatic cancer, with promising results

Researchers discover a potential therapeutic target for deadly cancer

Date: April 17, 2019

Source: Salk Institute

Summary: Advanced pancreatic cancer is often symptomless, leading to late diagnosis only after metastases have spread throughout the body. Now, researchers have uncovered the role of a signaling protein, called LIF, that may be the Achilles' heel of pancreatic cancer.

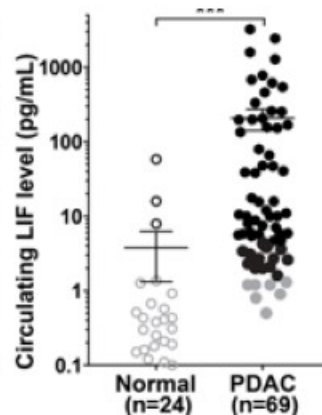


Credit: Salk Institute

Simoa LIF assay is more sensitive than Luminex

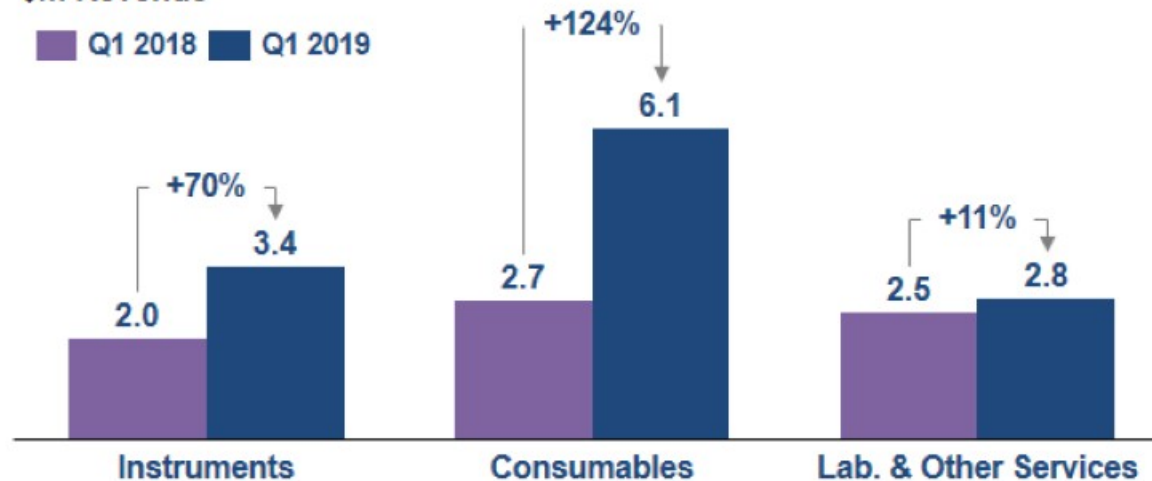
| | Luminex | | | | Simoa | | | |
|-------|------------|-------|-------|---------|------------|---------|---------|----------|
| | LIF(pg/mL) | FI-1 | FI-2 | Avg. FI | LIF(pg/mL) | AEB-1 | AEB-2 | Avg. AEB |
| Blank | 0 | 64 | 70 | 67 | 0 | 0.0143 | 0.0133 | 0.0138 |
| S1 | 6 | 87 | 78 | 82 | 0.11 | 0.0252 | 0.0200 | 0.0226 |
| S2 | 24 | 141 | 132 | 137 | 0.44 | 0.0420 | 0.0403 | 0.0411 |
| S3 | 98 | 323 | 303 | 313 | 1.76 | 0.1017 | 0.1121 | 0.1069 |
| S4 | 391 | 914 | 903 | 908 | 7.03 | 0.3762 | 0.3867 | 0.3815 |
| S5 | 1563 | 2745 | 2802 | 2773 | 28.13 | 1.5377 | 1.4952 | 1.5164 |
| S6 | 6250 | 7260 | 7129 | 7194 | 112.50 | 5.4800 | 5.6538 | 5.5669 |
| S7 | 25000 | 12974 | 12668 | 12821 | 450.00 | 21.3583 | 20.4018 | 20.8801 |

LOD: 0.048 pg/mL

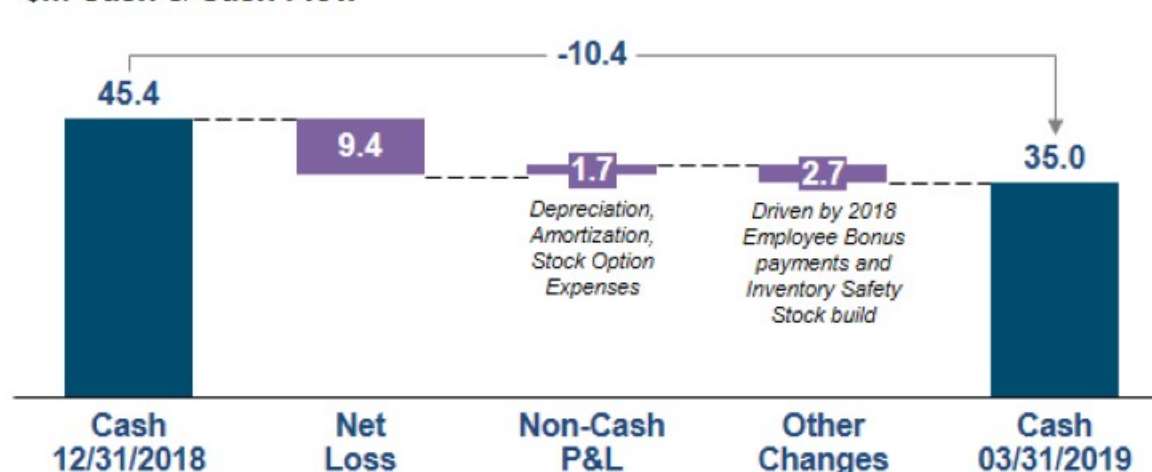


| in \$m | Actuals | Prior Year | |
|-----------------------------|--------------|-------------|-------------|
| | Q1 2019 | Q1 2018 | Var. |
| Product Revenue | 9.5 | 4.7 | +4.8 |
| <i>Growth vs. PYR</i> | +101% | | |
| Lab/Services | 2.8 | 2.5 | +0.3 |
| <i>Growth vs. PYR</i> | +11% | | |
| Collaboration | - | 0.3 | -0.3 |
| Total Revenue | 12.3 | 7.5 | +4.8 |
| <i>Growth vs. PYR</i> | +64% | | |
| Gross Profit | 6.0 | 3.2 | +2.8 |
| <i>Gross Margin %</i> | 48.7% | 42.2% | +650 bps |
| Operating Expenses | 15.4 | 10.3 | +5.0 |
| Loss from operations | -9.4 | -7.2 | -2.2 |

\$m Revenue



\$m Cash & Cash Flow



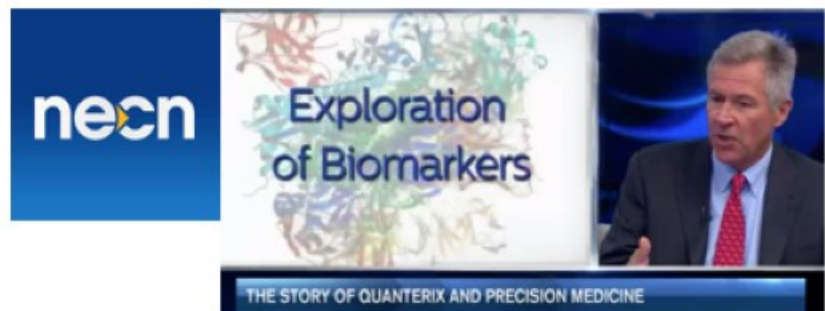
Poised to Disrupt Healthcare and Create Significant Value



Quanterix continued to build awareness for its mission to advance the science of precision health in mainstream press and trade media.

PPH 2018 - Amsterdam

MEDIA HIGHLIGHTS



The Story of Quanterix and Precision Medicine

Hrusovsky believes that in order to succeed, a business must change the world in a positive way. As he said, "We believe that by powering precision health, we are helping to eradicate the causes of Alzheimer's, ALS, and Parkinson's disease."

Forbes

ABC – Feb 2019

Bloomberg



Alzheimer's Might be Detectable With Biomarker Test 16 Years Before Onset of Symptoms

BOSTON BUSINESS JOURNAL



Bloomberg Radio Feb 2019



New technology could detect cancer in blood before symptoms develop

HUFFPOST



Neurofilament Light Chain Levels in Blood of Value as Biomarker of MS Activity and Treatment Response, Study Finds