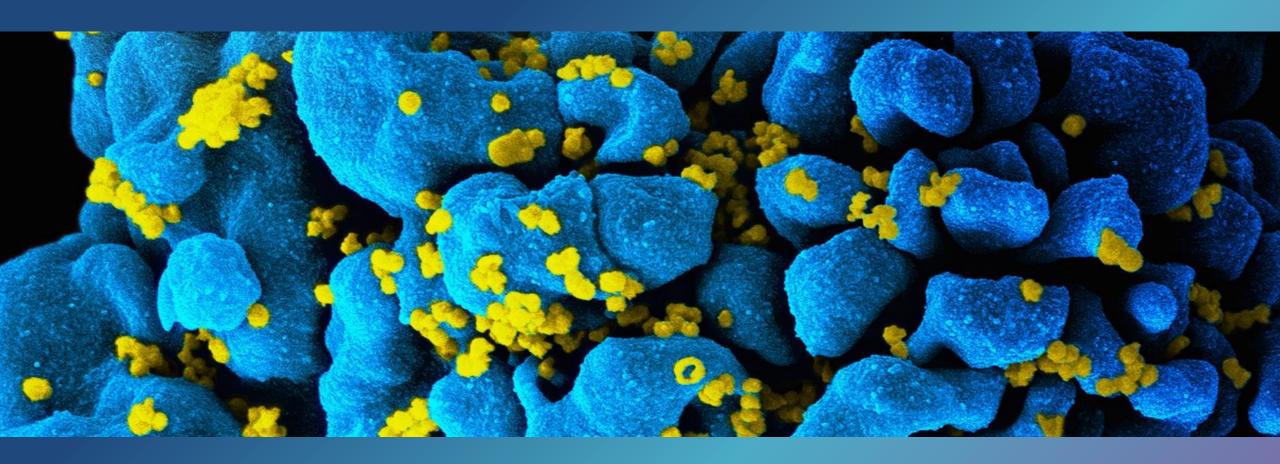
# Quanterix® The Science of Precision Health



BIO 2019 • June 5<sup>th</sup> 2019 Kevin Hrusovsky, Chairman & CEO

## Forward-Looking Statements



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.

# Today's Agenda



- 1. Compelling Strategic and Financial Opportunity
- 2. Compelling Milestones Achieved with Attractive Growth Catalysts
  - a. Record 2018 and Q1'19 Growth, Gross Margin and Growth Acceleration
  - b. \$50M Raise, New Head Quarters, China Launch, Consumable Expansion
  - c. SP-X Launch for Oncology and HD-X Launch Announced for Neurology
- 3. Revolutionizing Drug Development and Healthcare with Precision Health



# Today: Invasive and Late









# Tomorrow: Non-invasive and Early



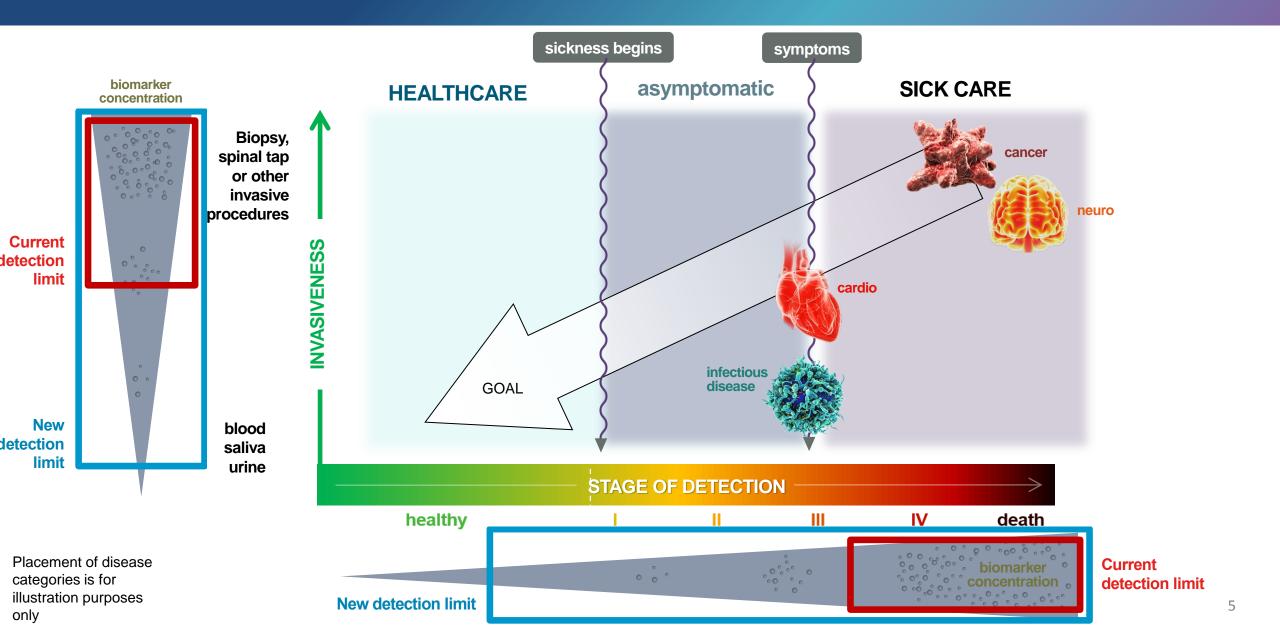






## Simoa Sees Health to Disease Continuum

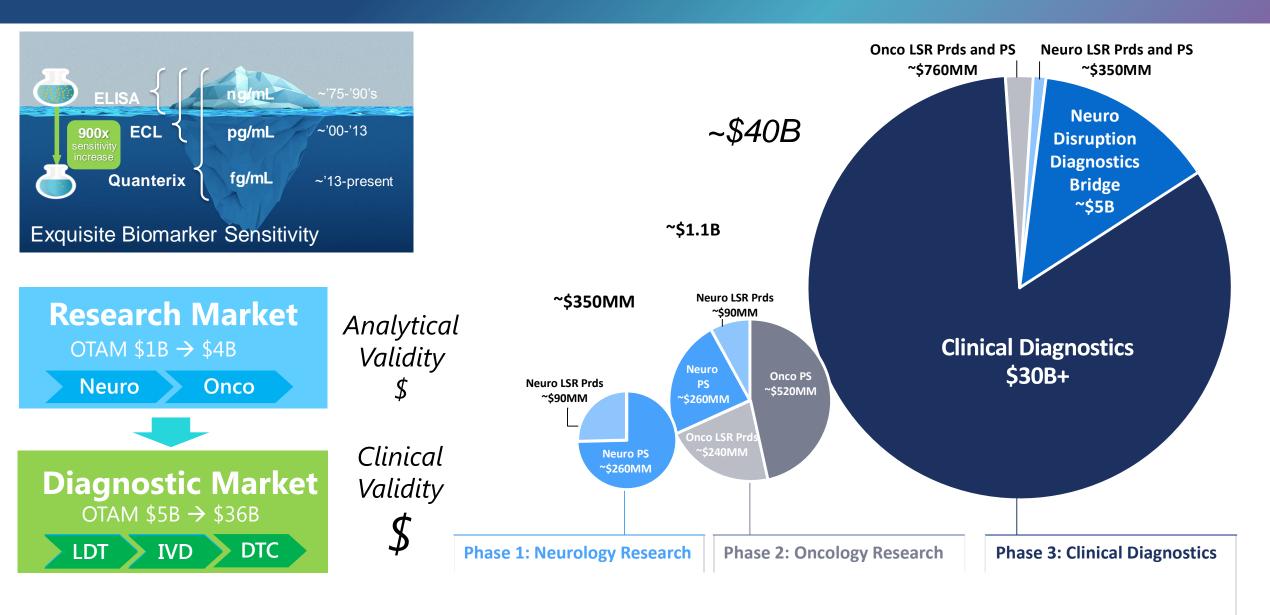




# Strategic Direction Remains Unchanged . . .

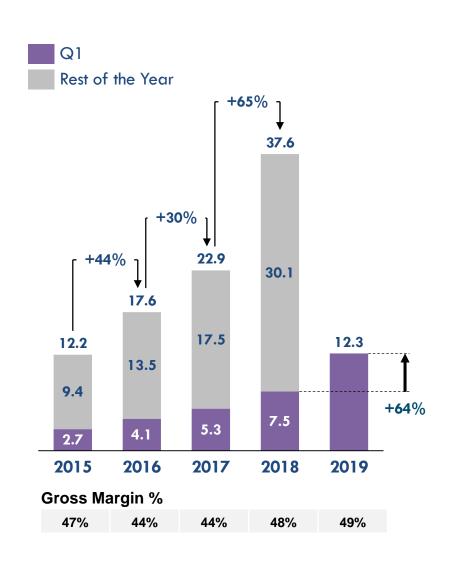
Capture major market opportunity through our disruptive technology





# Q1 2019 Growth Led by Consumables and Instruments



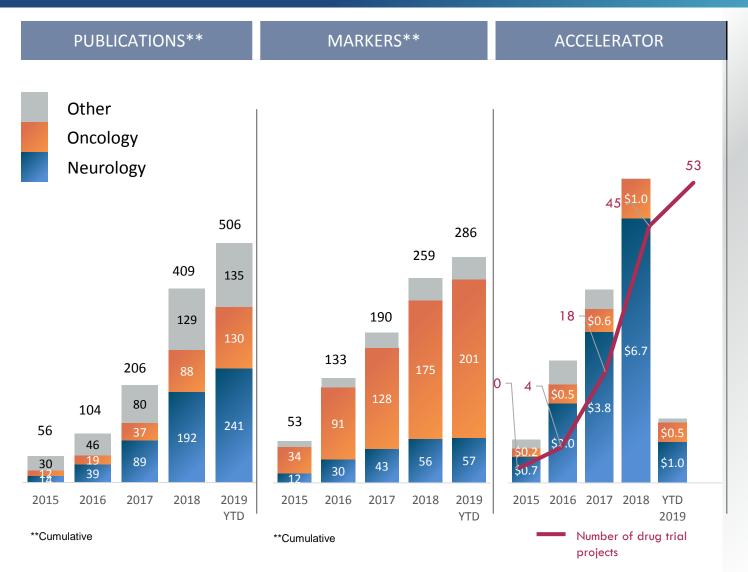


- 64% Growth, Total GM 49%, + 650bp
- High margin mix accelerating
- Consumables growth +124%



## Scientific Research is Driving Brand Awareness, Performance and Utilization



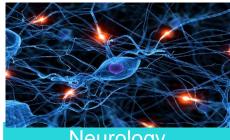




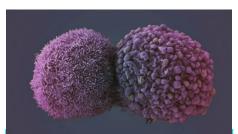
## Objectives 2019

### Large strides in Q1 towards securing our 2019 objectives





Neurology <10% Penetrated



Oncology 3x Neuro



Strategy
Dx via Biopharma



Financials



**Technology** 

High double digit growth with high utilization

Add 25 assays and globalize

Launch penetration in Oncology market

Immuno therapy focus

LDT/IVD partnerships

50 phase I, II, III trials

M&A: Ab's, Clinical Lab

LT Growth: 40%

Gross Margin: +300 bps

Instrument Growth: +25%

100x sensitivity by '21 YE

New frontier of medicine

Protein Translational Modifications

• +70% growth

- Launched SPX and 10 Simoa panels in Q2'19
- Impactful webinar, high utilization, early launch success
- Clinically validate Nf-L for 2<sup>nd</sup> ary endpoint & DP
- Partnerships progressing
- · Abbott blood screening

- Revenue growth +64%
- Gross Margin +650 bps
- Instrument growth +70%
- 40X defined
- Prototype developed

## **Quanterix Product Offerings**



Instruments	Assay kits	Services
-------------	------------	----------









HD-1 / HD-X Q4'19

SR-X

SP-X Q2'19

Plate

Bead

Accelerator

Floor-standing integrated

Assay prep and detection

400+ publications

Benchtop semiautomated assay prep Simoa planar assay

Benchtop semiautomated assay prep

Multiplex capabilities

250+ assays

Homebrew kits

Singleplex and multiplex \

Contract research & testing

Custom assay development & reagent production

**CLIA** and **LDT** capabilities

# We are Addressing a Significant Unmet Need in Drug Development FDA Announces Office of Drug Evaluation Science - ODES

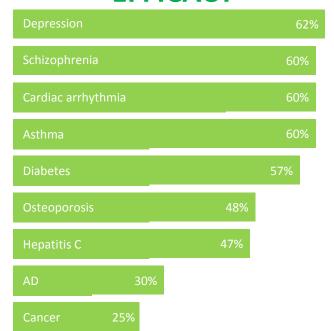


#### DRUG PERFORMANCE

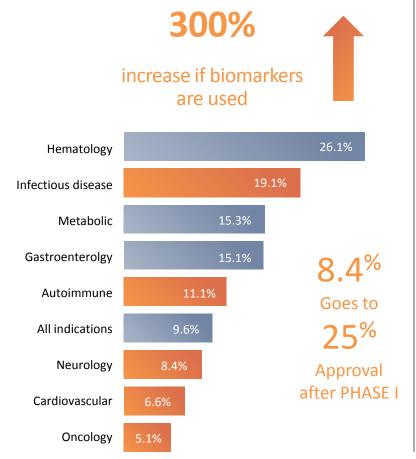
### **TOXICITY**

Adverse drug events are a substantial cause of Death in USA

### **EFFICACY**

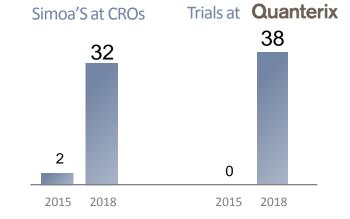


#### PROBABILITY OF DRUG APPROVAL

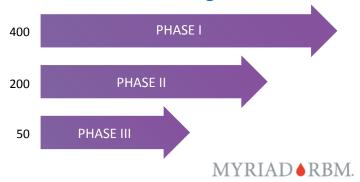


#### Probability of phase III approval after Phase 1 approval

#### VALIDATION OF SIMOA IMPACT



# 650 clinical trials with Simoa at single CRO



## Blue Chip Customers

· Cimi





# Other









✓ BioChek













































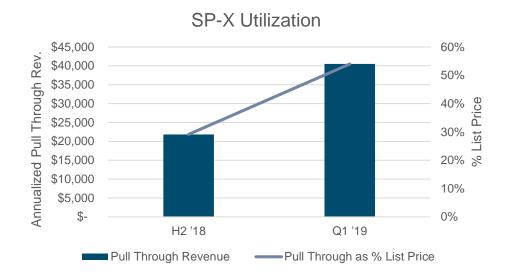


## SP-X Launch in Oncology

## **Exceptional Utilization at Key Accounts**







- **5** SP-Xs sold, **+86%** growth in utilization for top 3 customers
- 10 new multiplex panels across **52** immune & oncology biomarkers
- Equivalent sensitivity with HD-1

#### Robust Demo Program

























U NOVARTIS













# Launching Simoa HD-X Analyzer ahead of schedule Delivering digital biomarker disruption that you can count on





Re-engineering ultra-sensitivity, leveraging years of experience

Major productivity improvements, greater flexibility and unparalleled sensitivity

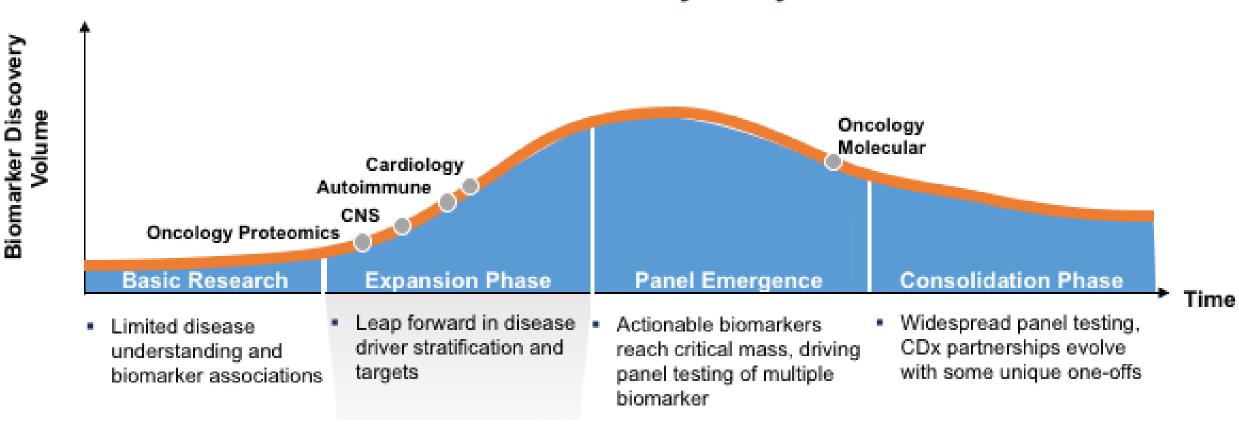
**Best-in-class assay performance** across a **broad assay menu** to empower biomarker research and accelerate drug development.

**Regulatory Compliance** – Enables 21CFR Part 11 compliance with streamlined run reports and user management

- Taking orders, shipping Q4
- Trade-in program available



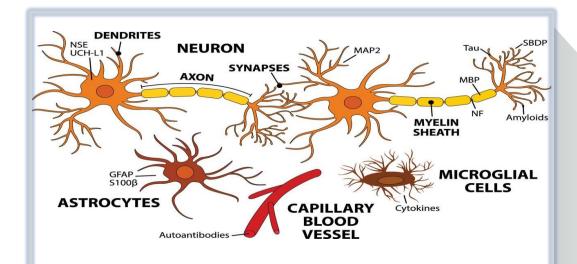
# **Biomarker Discovery Lifecycle**



### **Unmet Need**

### Development of neurological disease diagnostics has been slow due to disease complexity.





Neuro health has become national Health Priority due to veteran PTSD, opioid addictions, AD demographics mental health issues, and healthcare burden of neurological conditions Biopharma industry discouraged by lack of returns and use of subjective cognitive endpoints

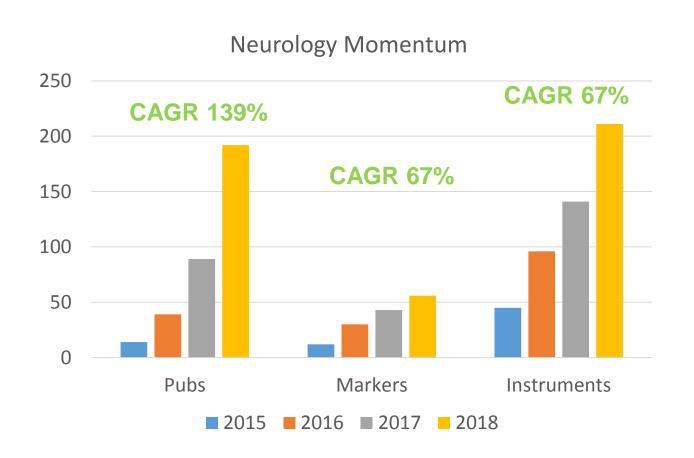


FDA issuing guidance to enable use of biomarkers in drug trials for early stage disease cohorts

Biopharma deploying biomarker approaches to trial design

## **Neurology Momentum Fueling Growth**





- Nf-L showing tremendous potential for key clinical applications in top peer-reviewed pubs; FDA interest
- Better AUC measuring AB40/42 than breakthrough device exemption for Alzheimer's
- Promise of blood-based neurology tests based on publications (CNN, Forbes, Bloomberg, Washington Post, GMA, etc.)
- Deep adoption by leading academic and pharma with strong network of world-leading KOLs















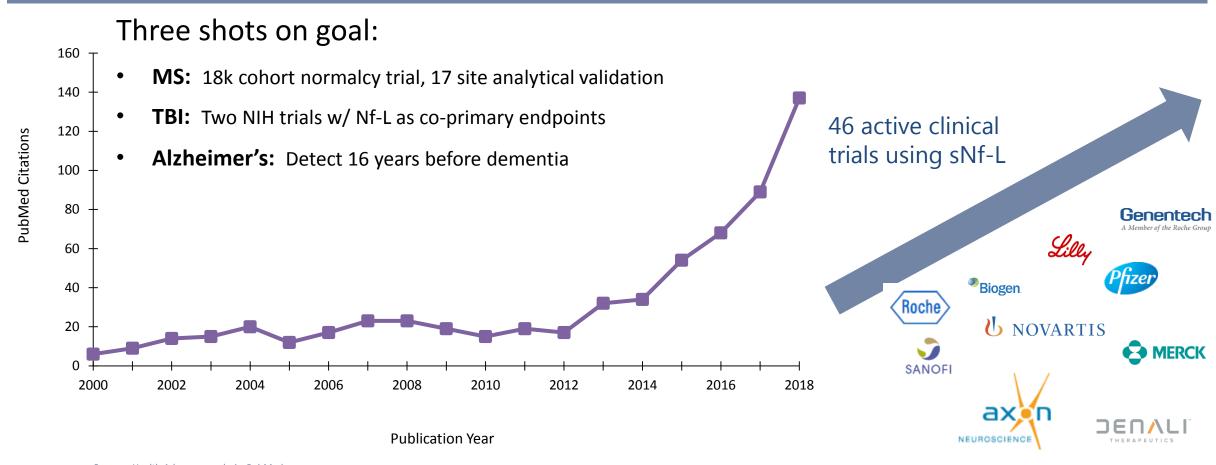




## NF-L Rapidly Expanding as Best In Class Neuro Biomarker



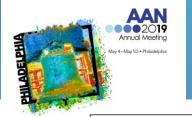
#### RESEARCH PUBLICATIONS ON NFL



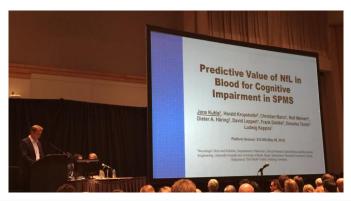
Source: Health Advances analysis, PubMed.

## AAN 2019: Explosive momentum of Simoa NfL

Search





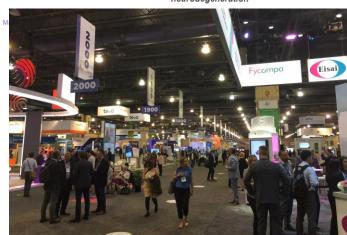


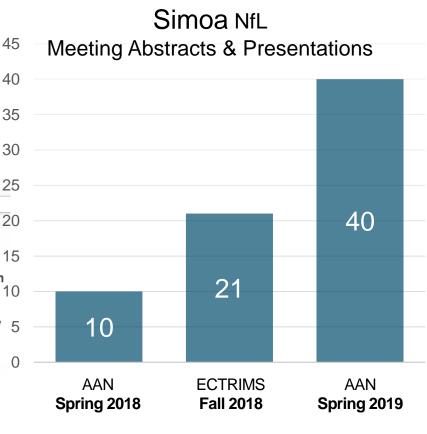


HOME SERVICES NEWS EDUCATION ABOUT US

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 (Nf-L) as a potential diagnostic and prognostic biomarker for





5 Phase III trials across >1600 MS patients demonstrate clinical utility of Simoa NfL

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

Peter A. Calabresi, <sup>1</sup> Jens Kuhle, <sup>2</sup> Douglas L. Arnold, <sup>5</sup> R. Philip Kinkel, <sup>4</sup> Ludwig Kappos, <sup>2</sup> Carol M. Singh, <sup>5</sup> Dipen Sangurdekar, <sup>5</sup> Carl De Moor, <sup>5</sup> Bob Engle, <sup>5</sup> Ray Su, <sup>5</sup> Aaron Deykin, <sup>5</sup> Elizabeth Fisher, <sup>5</sup> Alfred Sandrock, <sup>5</sup> Bernd C. Kleseler, <sup>6</sup> Richard A. Rudick, <sup>5</sup> Tatlana Plavina<sup>6</sup>

Department of Neurology, Johns Hopkins University School of Medicine, Baltimore, MD, USA: \*Neurologic Clinic and Policilinic, Departments of Medicine, Biomedicine and Clinical Research, University Hospital Basel, Basel, Switzerland: \*Montreal Neurological Institute, McGill University, Montreal, CC, Canada: \*Department of Neurosciences, University of California, San Diego, CA, USA: \*Biogen 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<sup>1</sup>, Finn Sellebjerg<sup>2</sup>, Hans-Peter Hartung<sup>3</sup>, Douglas Arnold<sup>3</sup>, Mark S. Freedman<sup>3</sup>, Douglas Jeffery<sup>3</sup>, Aaron Miller<sup>7</sup>, Keith R. Edwards<sup>3</sup>, Carol M. Singh<sup>3</sup>, Ih Chang<sup>3</sup>, Zhang Ren<sup>3</sup>, Dipen Sangurdekar<sup>7</sup>, Bing Zhu<sup>3</sup>, Devangi Mehta<sup>3</sup>, Pei-Ran Ho<sup>3</sup>, Nolan Campbell<sup>9</sup>, Michael Edwards<sup>9</sup>, Elizabeth Fisher<sup>9</sup>, Bernd C. Kieseier<sup>9</sup>, Richard A. Rudick<sup>9</sup>, Tatiana Playina<sup>3</sup>

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

Jeffrey Cohen<sup>1</sup>, Ludwig Kappos<sup>2</sup>, Nadia Tenenbaum<sup>3</sup>, Jackie Han<sup>3</sup>, Harald Kropshofer<sup>4</sup>, Davorka Tomic<sup>4</sup>, Je

<sup>1</sup>Cleveland Clinic, <sup>3</sup>Neurologic Clinic and Policlinic, Department of Medicine, Clinical Research, Biomedicine and Biomedical Engineering, University Hospital and University of Basel, <sup>3</sup>Novartis Pharmaceuticals Corporation, <sup>5</sup>Novartis 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.

This post hoc analysis was based on data from patients who received fingolimod 0.5 mg once daily or placebofinetrefor 0.1-a (FN) 30 µg once weekly in pivotal studies (24-month FREEDORIV2-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 NIL assessments at baseline, end of core (EoC) 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 placeboffR9 group to fingolimod in the LONGTERMS. Pit. was measured using Single Molecule Array (SIMOA\*\*) immunoassay. The geometric mean change in NIL levels from baseline to EoS was analyzed using Wilcoxon signed-rank test.

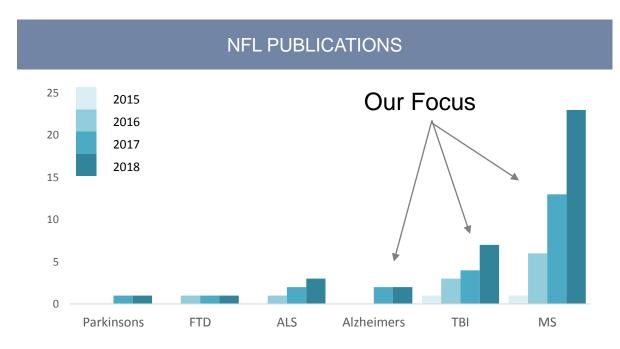
#### Results:

The mean exposure to fingolimod was 3483 days in the continuous group and 2822 days in the switched group. In the continuous group, baseline NIL levels of 33 pg/ml, were significantly reduced by approximately 40% at both EoC and EoS (20 pg/ml.; P-c0.0001 and P-0.0002, respectively). In the switched group, baseline NIL levels of 29 pg/ml. were reduced by 15% at EoC (25 pg/ml., P-o.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 RBMS.

## Emerging Clinical Biomarker: Neurofilament Light (NfL)



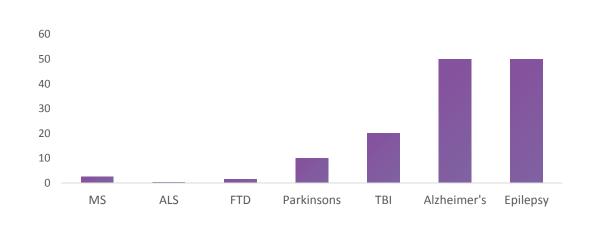


#### STUDIES CONFIRM NFL CLINICAL UTILITY:

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

Majority of published data obtained with Simoa NfL





#### **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

# Poised to Disrupt Healthcare and Create Significant Value



	Differentiator	Value
Category-defining; Unrivaled Sensitivity / Technology	Best in Class	Disrupt
market  Methodical market penetration strategy to reward investors	\$3B to \$38B	New Answers
3 DNA – RNA - Protein; Better linked to Disease / Health		Holy Grail
Quanterix 4 Validation: 19/20 top pharma, PPH, 800+ trials	500+ pubs All Areas	Proven
execution  Growth & Value; Razor – razor blade, \$150M invested	Product Launches	Rapid Growth
6 Low Risk / Solid Return + Uber Return Prospect		Retail
7 Track Record for Commercializing Disruption		Lynchpin