

J.P. Morgan Healthcare Conference 2021

Kevin Hrusovsky, Chairman, CEO & President January 13, 2021

Forward-Looking Statements & Non-GAAP Financial Measures

This presentation contains forward-looking statements within the meaning of the Private Securities Litigation Reform Act of 1995. Words such as "may," "will," "expect," "plan," "anticipate," "estimate," "intend" and similar expressions (as well as other words or expressions referencing future events, conditions or circumstances) are intended to identify forward-looking statements. These forward-looking statements include the Company's estimated Q4 and full year 2020 financial performance, which is preliminary and unaudited and is subject to completion of the Company's normal guarter and year-end close procedures. These procedures and the audit of the Company's financial statements for the year ended December 31, 2020 are ongoing and could result in changes to such estimated information. Forward-looking statements in this news release are based on Quanterix' expectations and assumptions as of the date of this press release. Each of these forward-looking statements involves risks and uncertainties. Factors that may cause Quanterix' actual results to differ from those expressed or implied in the forward-looking statements in this press release are discussed in Quanterix' filings with the U.S. Securities and Exchange Commission, including the "Risk Factors" sections contained therein. Except as required by law, Quanterix assumes no obligation to update any forwardlooking statements contained herein to reflect any change in expectations, even as new information becomes available.

To supplement the Company's financial statements presented on a GAAP basis, the Company has provided certain non-GAAP financial measures. Management uses these non-GAAP measures to evaluate the Company's operating performance in a manner that allows for meaningful period-to-period comparison and analysis of trends in its business. Management believes that such measures are important in comparing current results with prior period results, and are useful to investors and financial analysts in assessing the Company's operating performance. The non-GAAP financial information presented here should be considered in conjunction with, and not as a substitute for the financial information presented in accordance with GAAP. Investors are encouraged to review the reconciliation of these non-GAAP measures to their most directly comparable GAAP financial measures set forth in the appendix of this presentation.



Today's Agenda



Vision and Strategy for Disruptive Value

- Di5 Methodology
- Proteomics Leadership
- Research + Diagnostics
- Leapfrog to Payors



2020 Compelling **Progress**

- COVID PIVOT
- RUO-Dx Highway
- Enabling AD Drugs
- Ecosystem Catalyzed



Powering Precision Health

- Neuro
- COVID
- Immunology
- Disruption



Proteomics Disruption Catalyzing Precision Health Transformation

Compelling Thesis to Lead Proteomics into Precision Health Era

Execution

2 to 3x Value Creation

RESEARCH & DISCOVERY

\$1B to \$20B TAM



- Validation: Pubs, Trials, Pharma Adoption
- 2014-20 Rev: \$0 \$85m; 55% CAGR* Public
- Raised \$350m; Valuation \$30m to \$2.0 B
- **ST & LT Growth Catalysts**
 - Increase penetration Neuro, COVID, Immuno
 - Menu expansion & drug / population trials
 - Scale consumables growth New COO
 - Expand into upstream Discovery tools

* 2017-2020 3Yr CAGR

Aspiration

10 to 15x Value Creation

DIAGNOSTICS & HEALTH SCREENS

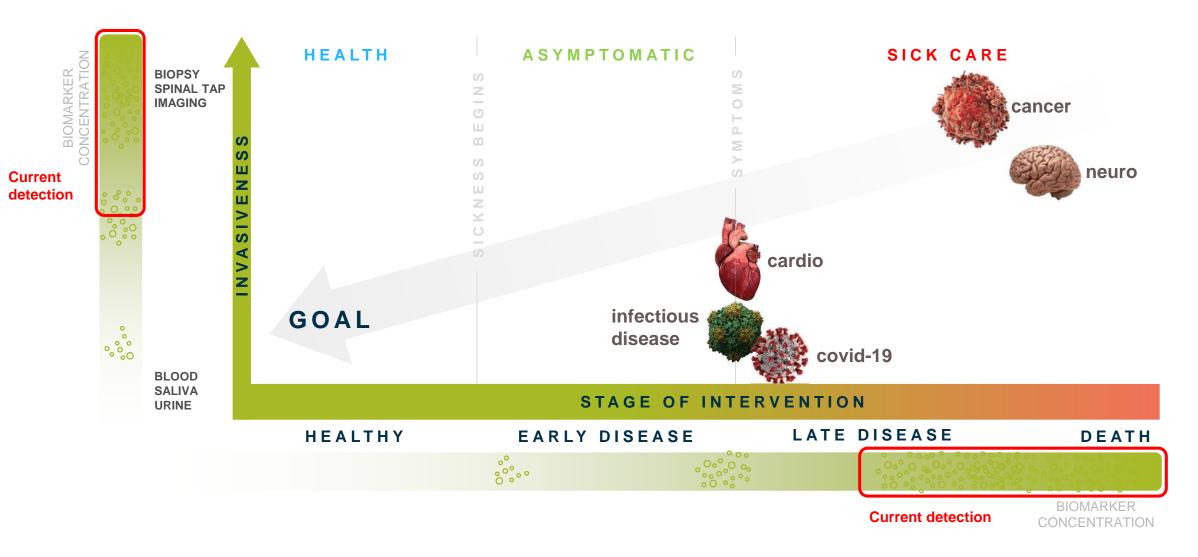
\$100B TAM



- Abbott, Siemens, Pharma Advances
- Payor Acceptance for Cost / Outcome
- NIH, RADx & FDA Ecosystem expansion (\checkmark)
- **ST & LT Growth Catalysts**
 - COVID EUA's
 - Neuro Tool Entry Leverage COVID
 - Liquid Biopsy Entry (Partner or Vertical)
 - Advance Payor "Outcome Leapfrog"

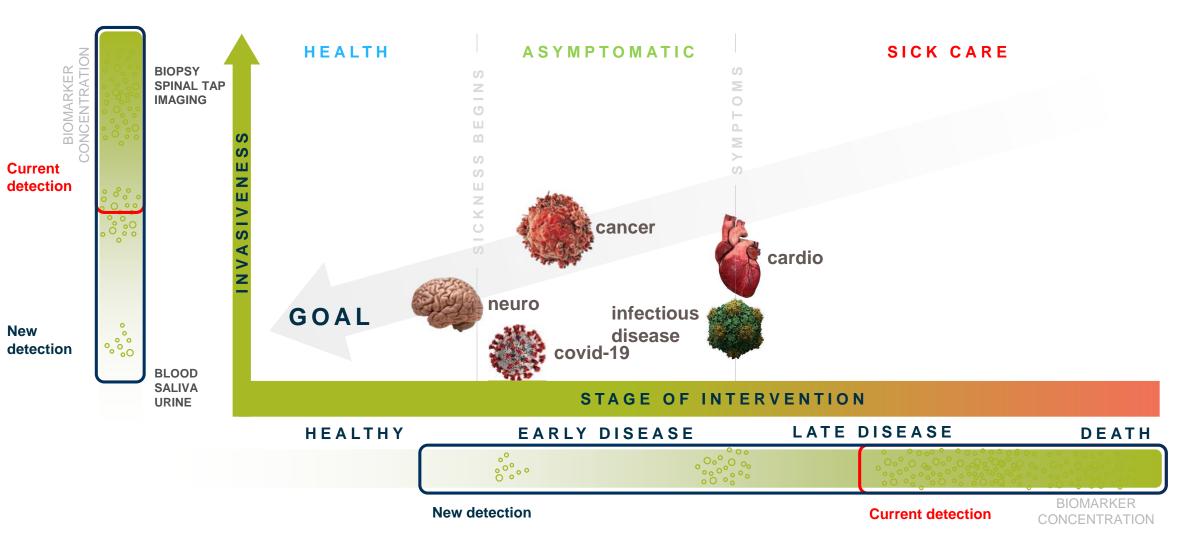
COVID **RADx**

Simoa® provides insight into Health to Disease Continuum



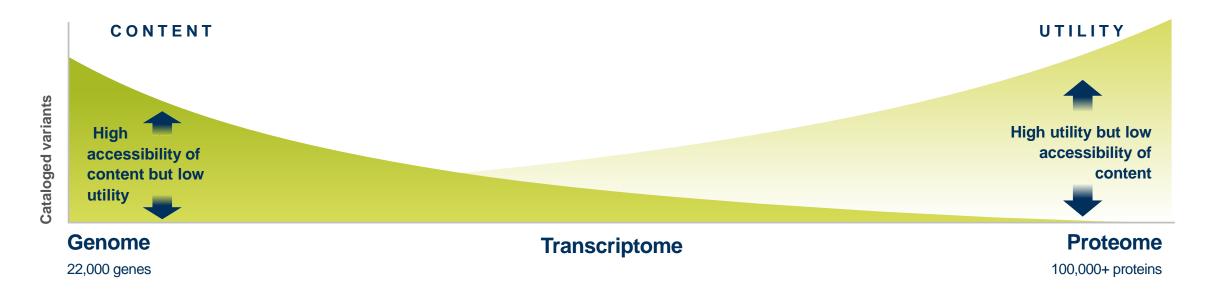


Simoa® provides insight into Health to Disease Continuum

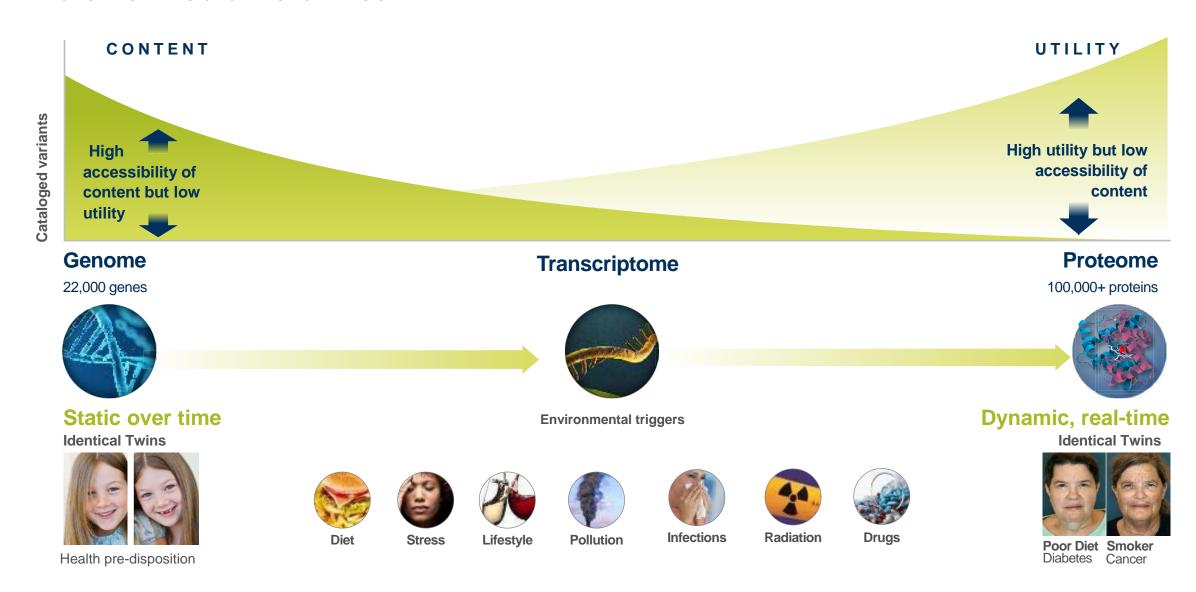




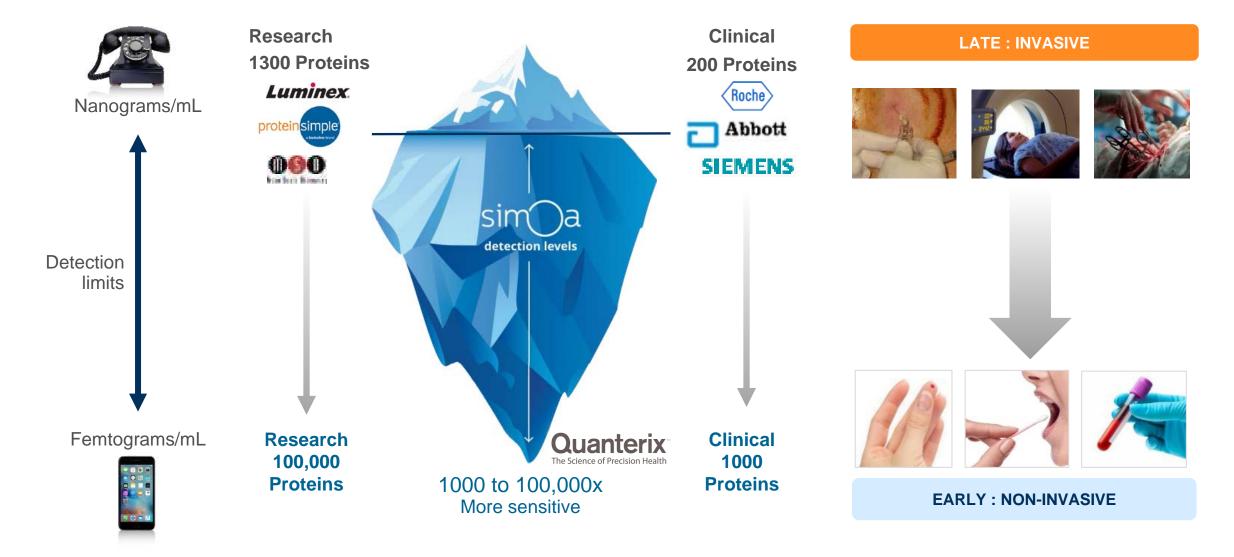
Proteins Predictive of Health



Proteins Predictive of Health



Digital Biomarkers Sensitivity Unlocking Proteomics



Disruption in sensitivity catalyzing TAM expansion & reallocation



Greater sensitivity enables...

- Less invasive, smaller samples
- New biomarkers & subtypes
- Multiplexing for disease specificity
- Single molecule to unravel heterogeneity

...which changes the healthcare paradigm:

- Early detection
- Health to disease continuum
- Homecare & frequent testing
- Disease progression monitoring
- Prevention, costs & outcomes

Revolution in Neuronal Biomarkers Measurement: CSF & Blood



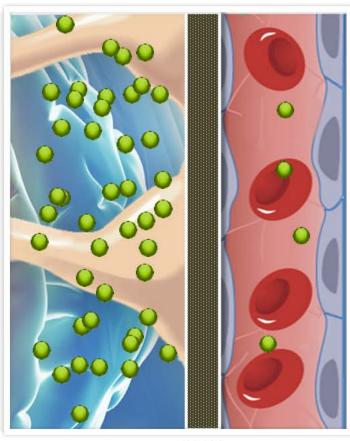
NfL

100 to 200 pg/ml

AB 42

451 pg/ml

Tau 87 pg/ml



Cerebral spinal fluid

BBB Blood vessels

Simoa Ultrasensitive detection required to quantify in blood

NfL

2 to 4 pg/ml

AB 42

Tau

4.70 pg/ml

1.65 pg/ml

Megatrends in Healthcare Catalyzing Proteomics for Precision Health



Rise of Home Care



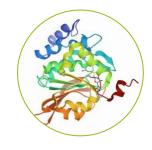
Employee, Student & Entertainment



Prevention & Asymptomatic



Inclusion & Equity



Proteomics Disruption





Decentralized Trials

2020 2021 2022 2023 2024 2025 2026 2027



Created Neurology Tools Foray into Immunology





2020

2021

2022

2023

2024

2025

2026

2027

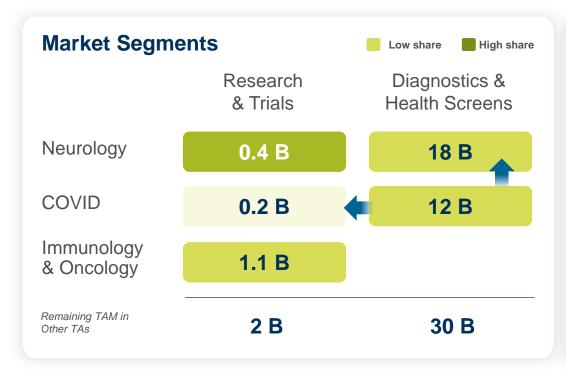




Created Neurology Tools Foray into Immunology

COVID Dx accelerates Neuro entry

Deepen Immunology research







2023 2024 2025 2026 2027 2020 2021 2022



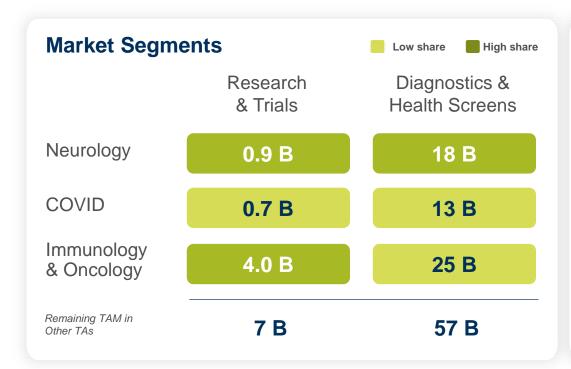
Post-COVID: **Accelerated Dx**

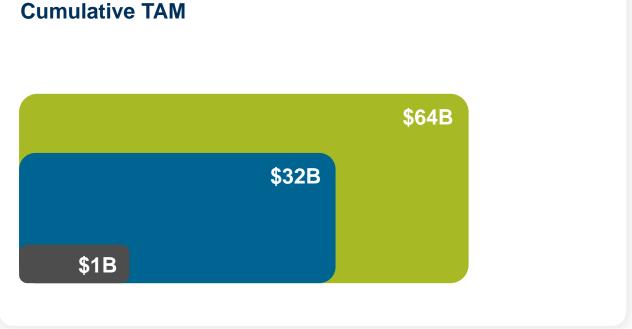
Created Neurology Tools Foray into Immunology

COVID Dx accelerates Neuro entry Deepen Immunology research



Proteomics liquid biopsy platform Trial decentralization & health screens





2025 2020 2022 2021 2023 2024 2026 2027





Created Neurology Tools Foray into Immunology

COVID Dx accelerates Neuro entry Deepen Immunology research

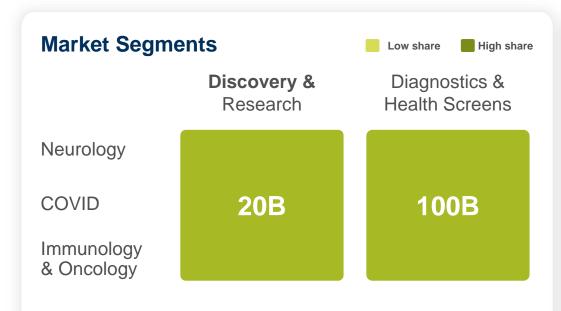


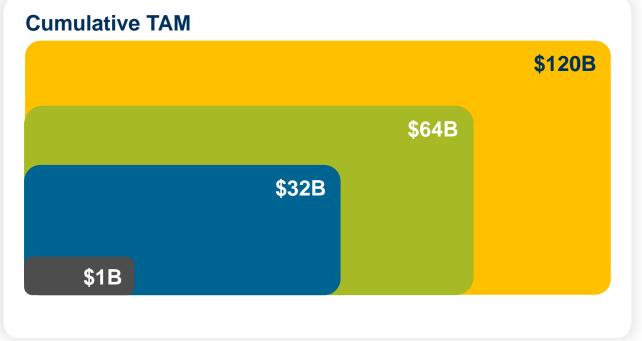
Disruptive growth: **PPH Megatrends**

Proteomics liquid biopsy platform Trial decentralization & health screens



Move upstream to catalyze biomarker discovery and increase protein flow





Near term priorities to catalyze & fortify proteomics leadership



COVID Vertical Fulfilment

Commercialize COVID EUA's

Drive Leading-Edge research including long-hauler research

Convert research into **Novel Differentiated Commercial Applications**



Neurology **Vertical**

Launch novel neuro multiplex; expand clinical trials and China

Continue Accelerator expansion and add **Neurology Drug rescue**

Launch Nf-L MS Minimum Residual Disease monitoring, Leverage Pharma data

Clinically validate **AD pre-symptomatic screening** and develop AD triage through pharma partnerships



Immunology Expansion

Launch new immunology & oncology biomarkers

Position SP-X as platform of choice for proteomics liquid biopsy



Disrupt Precision Health

Expand Payor Relationships for Health Screens – OUTCOME LEAPFROG Leverage Pharma and Payor relationships to expand DECENTRALIZED CLINICAL TRIALS

2020 Compelling Progress: Record Growth, Adoption & Value Ascent

COVID



- ✓ Serology & Antigen EUA's
- ✓ RADx \$20 million, NIH/FDA ties
- ✓ Payor Adoption Population studies
- ✓ Leading-edge research & capability

Neurology

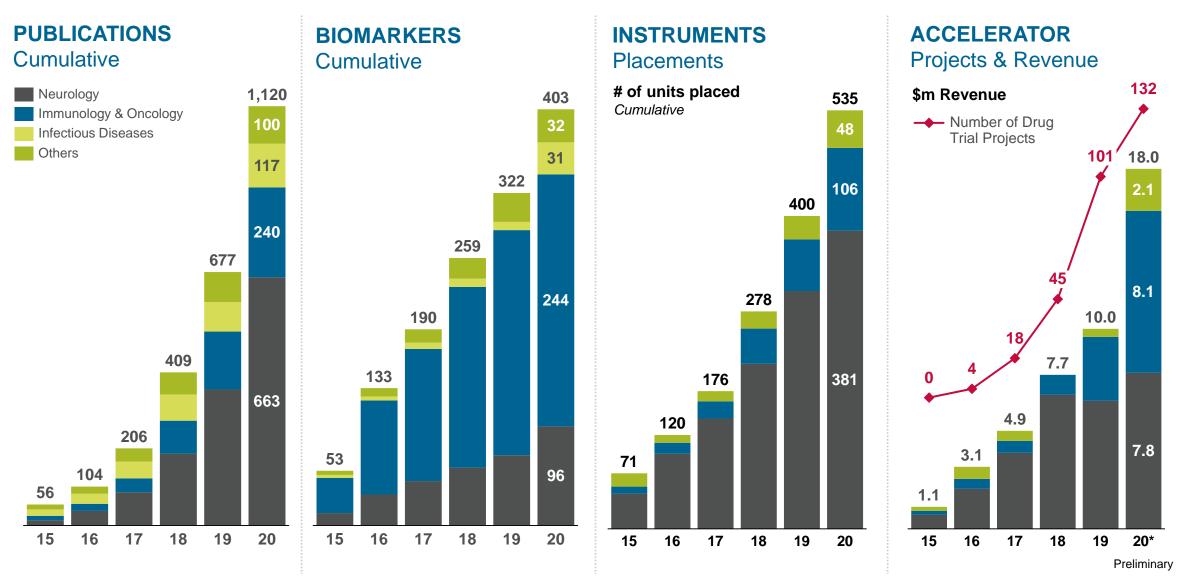


- ✓ Launched pTau-181 & Neuro panel
- ✓ Alzheimer's drug trials momentum
- MS Drug approvals with NfL
- √ ~48% HD-Fleet is HD-X & GROWTH



- Record publications & funding
- √ 100x Sensitivity pilot success
- ✓ Abbott License Agreement
- ✓ Lab Services Expansion & Growth

Scientific Validation driving Adoption





Superb Execution Continues

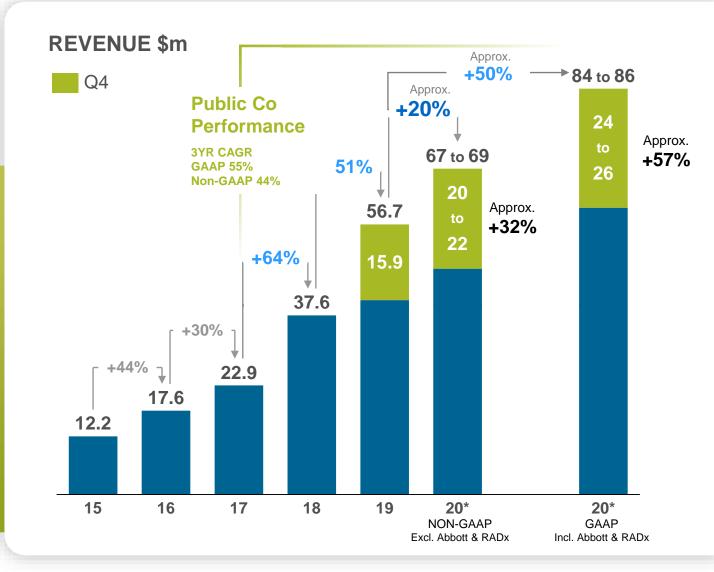
Strategic Growth Indicators

lncrease in Accelerator revenue 2019 to 2020, driven by immunology

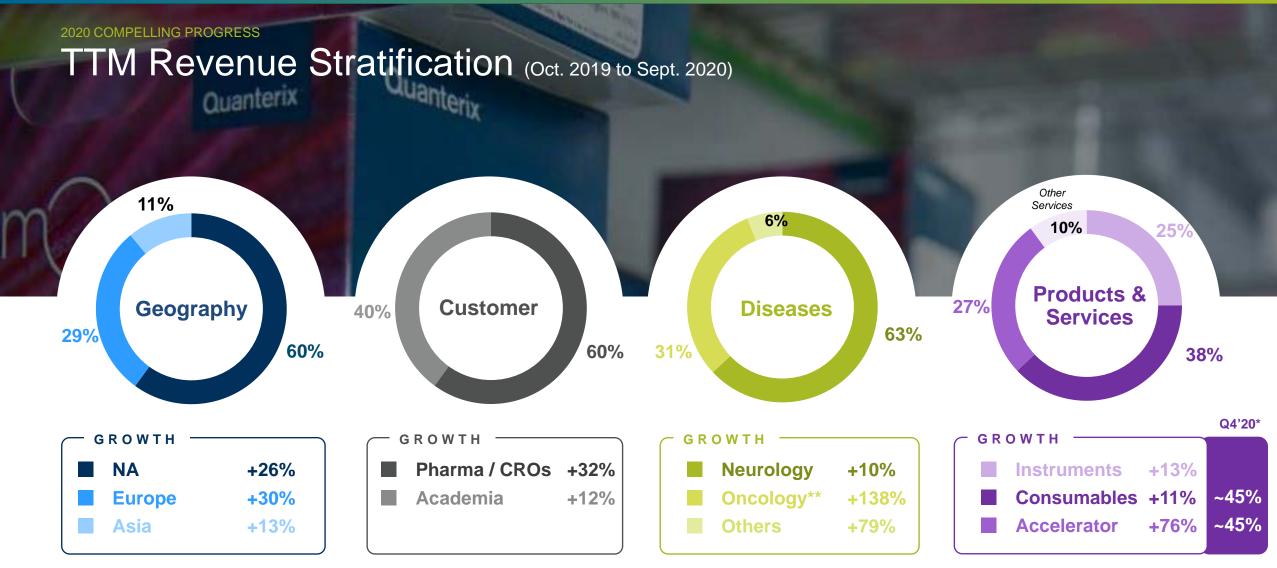
48% Proportion of HD-Fleet on new HD-X technology at YE2020

180 Instruments installed, with 93 HD-Xs (50%+ New) and 84 SR-Xs & SP-Xs

Consumables Revenue growth in Q4 2020 vs. Q4 2019 with utilization reverting to 55%+







Neurology back to 30-40% Growth in Q4*

All numbers exclude 1H FY19 \$1M 1x Study revenue



^{*} Q4 2020 and FY 2020 information are preliminary, unaudited estimates and are subject to the Company's normal quarter and year-end close procedures. These procedures and the audit of the Company's financial statements for the year ended December 31, 2020 are ongoing and could result in changes to such estimated information.

^{**} Incl. Immunology & Inflammation;

Delivered 2020 Objectives Plus, Despite Headwinds



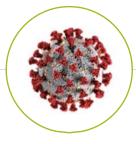
Neurology

Despite COVID headwinds . . .

Double-digit growth

48% HD-X IB at YE 2020

APAC Lab. Services partnership and expanded presence



COVID

Serology EUA

Antigen EUA

RADx Contract

Top-Payor Surveillance Studies



Strategy

Abbott License Agreement

AD Trials use of QTRX biomarkers

Several Diagnostic Accelerators across COVID and Neurology



Financials

FY Reported Revenue Growth ~50%

Non-GAAP Revenue* Growth ~20% despite COVID

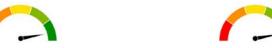
93 HD-X placements (50%+ New Installs)

84 New SR-X/SP-X placements



Technology

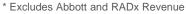
100x Sensitivity increase pathway defined and published











Objectives 2021

RUO 2019-2024 CAGR 30-40% - Diagnostics Pre-Mature and Incremental . . .

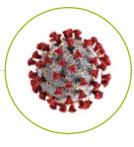


Neurology

Increases Biomarker penetration in Neuro Trials

65% HD-X IB at YE 2021

Measurable progress on Neuro Diagnostic Vertical & Payor – Pharma Drug Trials



COVID

Drive Antigen & Serology Commercialization

Deliver RADx Scale-Up

Support Leading-Edge COVID Research and develop novel solutions

Expand Payor Surveillance Studies



Immunology

Expand Assay Menu & accelerate publication flow

50 SP-X placements

Progress SP-X towards platform of choice for Liquid Biopsy LDTs

NIH – Payor – Pharma Longhauler & Drug Trials



Financials

RUO 5Yr CAGR 30-40% over 2019-2024

Diagnostics (COVID and Neuro) Pre-mature and Incremental

105 HD-X placements (60%+ New Installs)

140 New SR-X/SP-X placements



Platform

Operation Scale Quanterix (OSQ)

Define pathway to deploy 100x Sensitivity in the field

Expand Di5 PPH
Disruptive
Relationships

Our mission and vision to power precision health



1 Therapies



2 Prevention

Advocates - KOL's - Agencies - Investors



PATIENTS

We are on a mission to change the way in which healthcare is provided by giving researchers the ability to closely examine the continuum from health to disease.

In order to make this vision a reality, we formed a collaborative ecosystem by bringing together the most experienced management team, renowned scientists, industry leading investors and expert advisors, all united through the common goal of advancing the science of precision health.



Brain Health



COVID



Immunology and Oncology



Disruptive opportunities

Value Creation & Disruptive Growth Accelerators



Di5

1 Greater Purpose

FIVE STEPS TO
COMMERCIALIZING
DISRUPTIVE
INNOVATION

2 PPH Ecosystem

3 Culture & Leadership

Operation Scale Q

- PPH Inside
- Efficiency & Capacity
- Product launches
- Org & infrastructure

4 Strategy Roadmap

5 Superb Execution





Neurology poised for Value Creation Chain Reaction

Drug-Trials Penetration center-piece of near-term focus

Biomarker Discovery

Research led by academia and pharma



Currently underpenetrated Partnering with

Pharma to accelerate adoption (End-Points, Patient Stratification)















Imaging & Sample data support Clinical Validity, Utility and HECON

Access to Trial data accelerates and reduces cost for Dx Clinical Study

NfL pTau181 pTau217 Αβ42,Αβ40 tTau **GFAP**



UnitedHealth Group



Path to Dx further fuels Drug Development Biomarker Adoption in **Drug Development** Diagnostic Launch, Adoption & Reimbursement

Accelerators Driving Rapid Adoption for Transforming Drug Development

Research Institutions

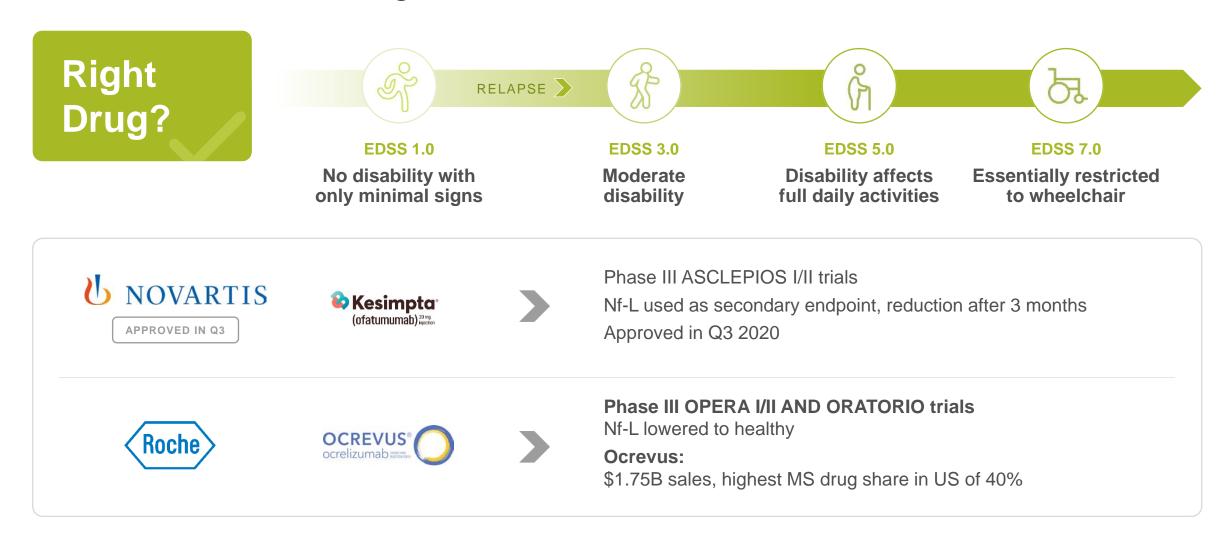


Biopharma Biogen Genentech Roche ODOVIE Scrum ADX & Bristol-Myers Squibb MERCK **∦**G∈n∈Works wren 冷泉港生物科技股份有限公司 Cold Spring Biotech Corp VIELABIO UNOVARTIS Achucarro MedImmune 2 Alnylam **B**IOMARIN DENAL chimera biotec Gen2 AMGEN GILEAD

CROs & Other



Serum Nf-L Powering Major MS Drug Trials \$22B Market – 16 drugs



AD Blood Biomarkers - Research to Clinical

Lilly & Quanterix Collaboration on Alzheimer's Disease Biomarkers

Research & Trials

< 0.5B TAM

Very early detection 16 yrs before dementia Blood p-Tau 217 / 181 / NfL Study: Blood test for Alzheimer's detects signs 20 years before memory, thinking falter Ken Alltucker USA TODAY Published 8:00 p.m. ET Ar. 28, 2020

0 9 2

A new blood test detected Alzheimer's disease as accurately as expensive brain seans or spinal taps, raising the possibility for a new, inexpensive option to

Researchers at the Alzheimer's Association International Conference on Tuesday presented the results of multiple studies of whether a blood test could distinguish Alzheimer's disease from other forms of dementia.

In one study published in JAMA, researchers said the blood test could could identify Alzheimer's disease and even detected signs of disease 20 years before

The New Hork Times **Bloomberg**





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

(CRN) — A simple blood test could predict if a publicant before symptoms begin, a new study finds.

fyr house of charges are a second of the protein could be an airly sign of the dissess second to the study outsides who has a least of the protein could be an airly sign of the dissess second to the study outsides blonday in the journal feature fedicine.

NS, is a "marker in the book whom glass an indicator of herve cellibros in the book," exp. more belief carrage you have," he said.







AD Blood Biomarkers - Research to Clinical

Lilly & Quanterix Collaboration on Alzheimer's Disease Biomarkers

Research & Trials

< 0.5B TAM

Very early detection 16 yrs before dementia

Blood p-Tau 217 / 181 / NfL

HEALTH

Study: Blood test for Alzheimer's detects signs 20 years before memory, thinking falter

Ken Alltucker USA TODAY Published 8:00 p.m. ET Ar. 28, 2020

0 💆 🖼

A new blood test detected Alzheimer's disease as accurately as expensive brain seass or spinal taps, raising the possibility for a new, inexpensive option to diagnose the most common form of dementia, researchers said.

Researchers at the Alzheimor's Association International Conference on Tuceday presented the results of multiple studies of whether a blood test could distinguish Abrheimer's disease from other forms of dementia.

In one study published in JAMA, researchers said the blood test could could identify Abrheimer's disease and even detected signs of disease 20 years before

The New Hork Times

Bloomberg











Diagnostics & Health Screens

~\$11B TAM

Relevance to Alzheimer's treatment

CSF p-Tau 181 modulated by aducanumab

HEALTH AND SCIENCE

Biogen's stock jumps 42% after FDA staff says it has enough data to support approving Alzheimer's drug





Hope for therapeutic with donanemab

BUSINESS | HEALTH CARE | HEALTH

Lilly Alzheimer's Drug Helped Patients in Small Trial

Donanemab slowed cognitive and functional decline of patients with mild Alzheimer's by 32% compared with placebo



AD Blood Biomarkers - Research to Clinical

Lilly & Quanterix Collaboration on Alzheimer's Disease Biomarkers

Research & Trials

< 0.5B TAM

Diagnostics & Health Screens ~\$11B TAM

Very early detection 16 yrs before dementia

Blood p-Tau 217 / 181 / NfL

WEALTH

Study: Blood test for Alzheimer's detects signs 20 years before memory, thinking falter

Ken Alltucker USA TODAY Published 8:00 p.m. ET Au. 28, 2020

O 💆 🖼

A new blood test detected Alzheimer's disease as accurately as expensive brain seans or spinal taps, raising the possibility for a new, inexpensive option to diagnose the most common form of dementia, researchers said.

Researchers at the Alzheimor's Association International Conference on Tuesday presented the results of multiple studies of whether a blood test could distinguish Alzheimer's disease from other forms of dementia.

In one study published in JAMA, researchers said the blood test could could identify Alzheimer's disease and even detected signs of disease 20 years before

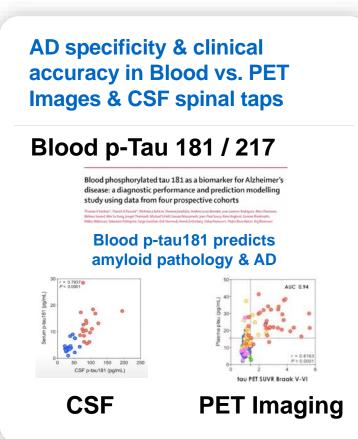
The New York Times

Bloomberg









Relevance to Alzheimer's treatment

CSF p-Tau 181 modulated by aducanumab

HEALTH AND SCIENCE

Biogen's stock jumps 42% after FDA staff says it has enough data to support approving Alzheimer's drug





Hope for therapeutic with donanemab

BUSINESS | HEALTH CARE | HEALTH

Lilly Alzheimer's Drug Helped Patients in Small Trial

Donanemab slowed cognitive and functional decline of patients with mild Alzheimer's by 32% compared with placebo



Recent excitement in liquid biopsy points to future potential of "liquid MRI"

Liquid biopsy for screening



JUN 2016

FDA approved first liquid biopsy test (lung cancer)

DEC 2019

(A) GUARDANT receives expanded Medicare coverage for solid tumor test

MAR 2020

* natera receives favorable Medicare reimbursement for organ transplant test

JUN 2020

INVITAE acquires ARCHER) in a \$1.4B deal

SEP 2020

illumina acquires GRALL at \$8.0B valuation

OCT 2020

SCIENCES acquires Thrive. at \$2.15B valuation

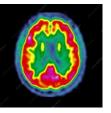
"Liquid MRI" for brain



Approval of first Alzheimer's Drug likely will be the tipping point for similar value acceleration and adoption momentum

Neurology Reimagined





CSF & PET Imaging

Αβ42	tTau	pTau181	
Nf-L	BDNF	SNAP-25	
α-synuclein	GDF-15	IL-6	
GFAP	Neurogranin	NPTX2	
S100b	sTREM2	TDP43	
VILP-1	YKL-40	MMP-3	
Αβ37/38/40	Osteopontin	IGFBP7	
	NSE	TRAIL	

Alzheimer's Parkinson's **Multiple Sclerosis Traumatic Brain Injury Lewy Body Dementia Huntington's** Stroke ALS FTD

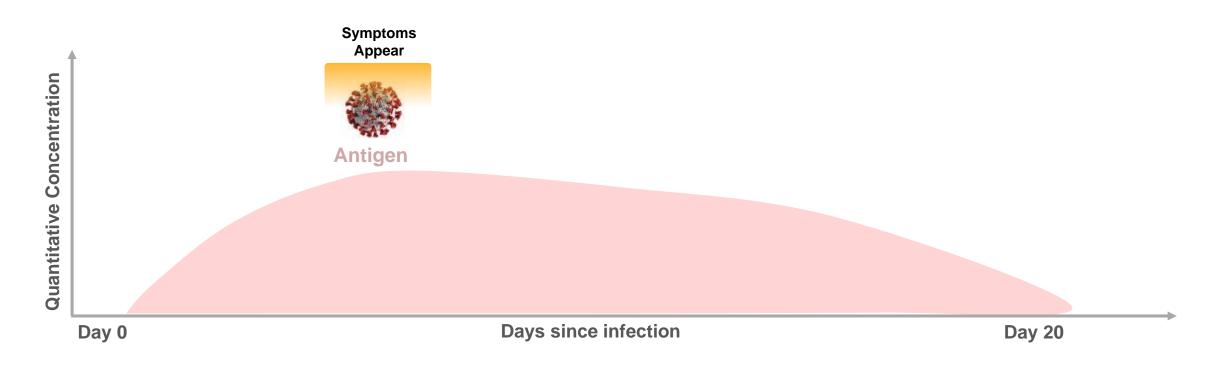


Blood-based SIMOA Neuro A	Assays
---------------------------	--------

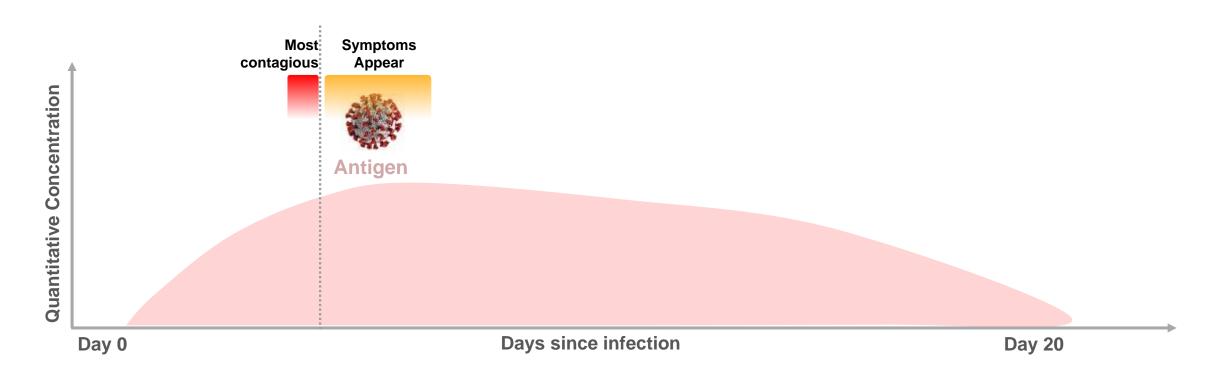
Αβ42,Αβ40	tTau	pTau181/217	SIEMENS
Nf-L	BDNF	SNAP-25	Roche
α-synuclein	GDF-15	IL-6	Abbott
GFAP	Neurogranin	NPTX2	BIOMÉRIEUX
S100b	sTREM2	TDP43	Biogen
VILP-1	YKL-40	MMP-3	b NOVARTIS
Αβ37/38	Osteopontin	IGFBP7	Lilly Eisai
	NSE	TRAIL	Roche

- Developed in Serum and/or Plasma Simoa catalog assay kits
- Developed in Serum and/or Plasma Simoa homebrew assays

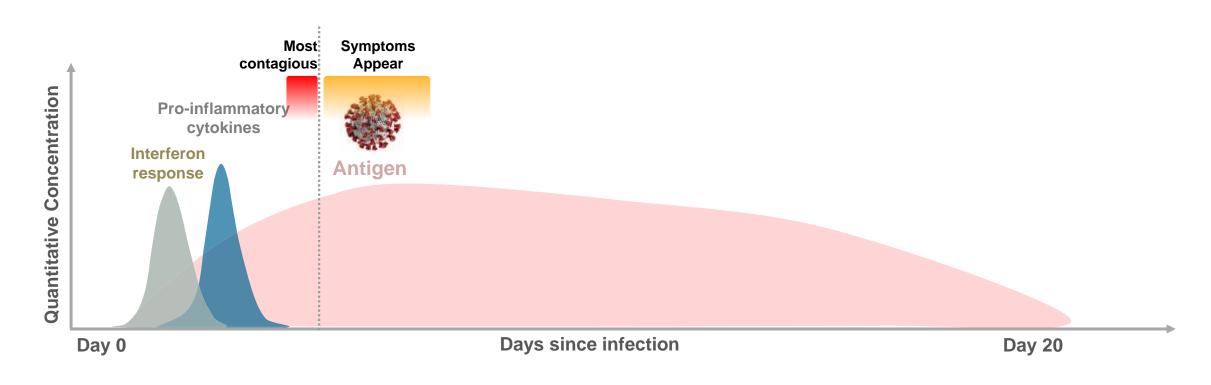
Roadmap to High-Definition picture of SARS-CoV-2 Infection



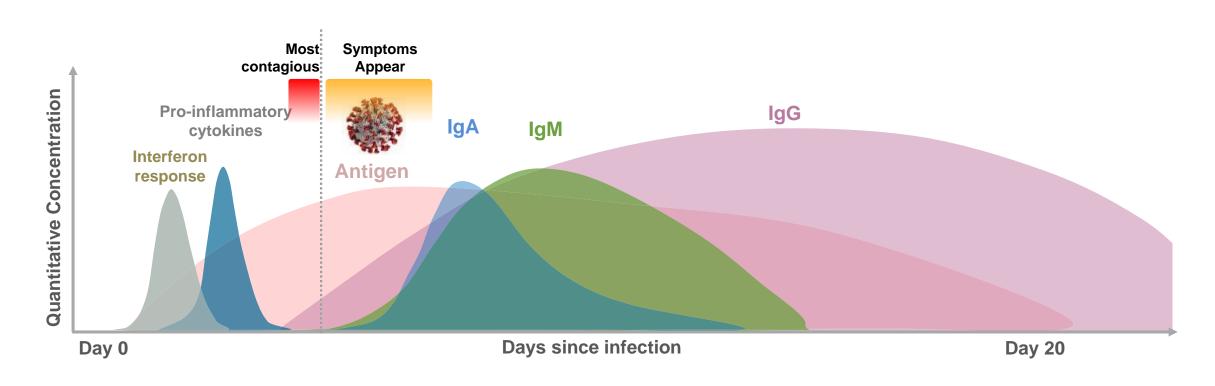






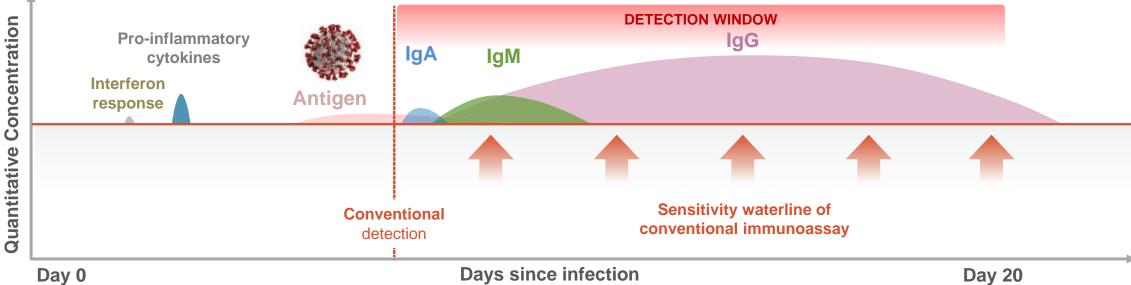




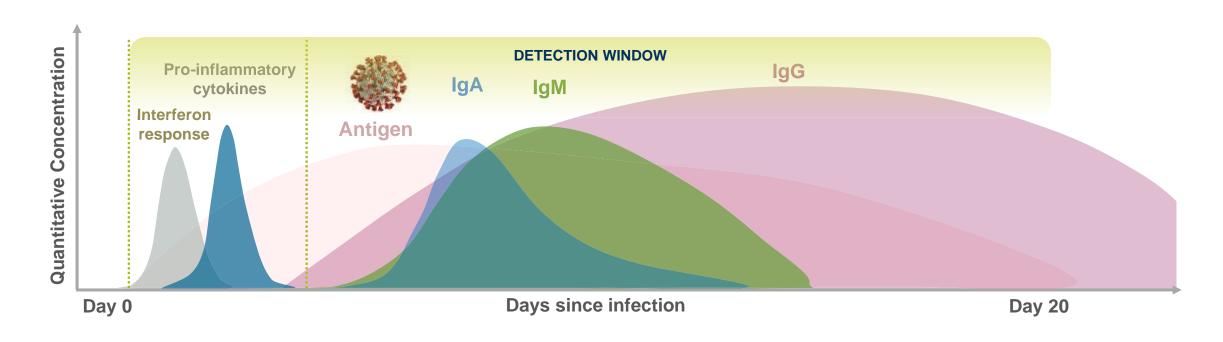


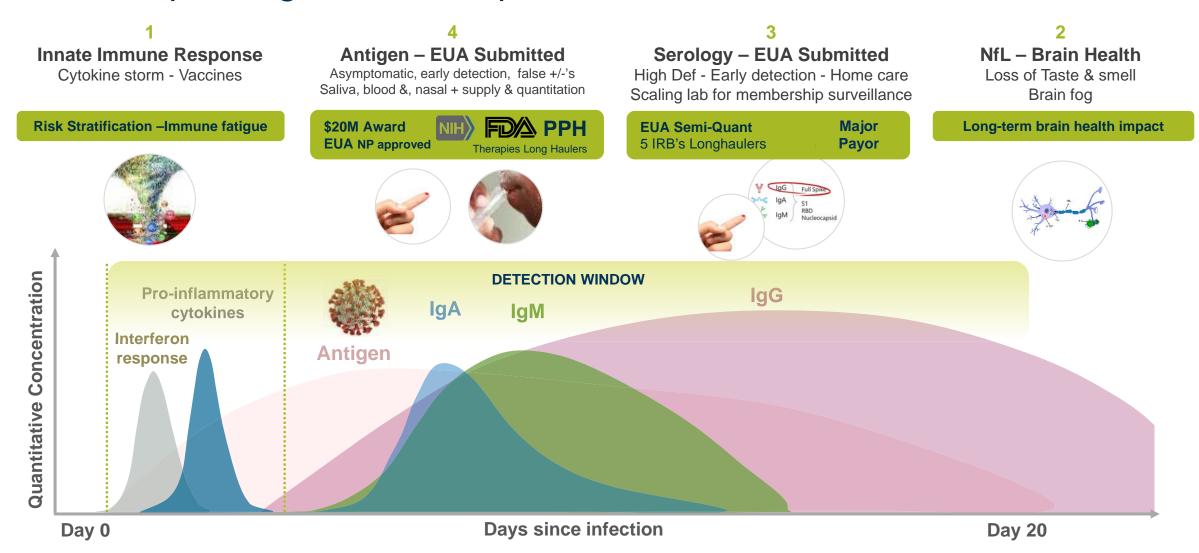






NEW TESTING PARADIGMDisruption Enabled By Ultra Sensitivity Digital ELISA





1

Innate Immune Response
Cytokine storm - Vaccines

Risk Stratification –Immune fatigue

4

Antigen – EUA Submitted

Asymptomatic, early detection, false +/-'s Saliva, blood &, nasal + supply & quantitation

\$20M Award EUA NP approved



3

Serology – EUA Submitted

High Def - Early detection - Home care Scaling lab for membership surveillance

EUA Semi-Quant 5 IRB's Longhaulers

Major Payor 2

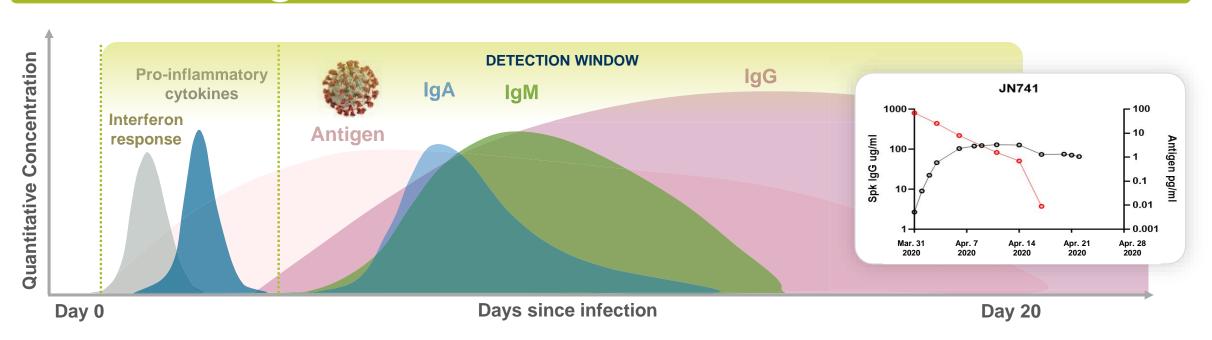
NfL – Brain Health Loss of Taste & smell

oss of Taste & smell.
Brain fog

Long-term brain health impact

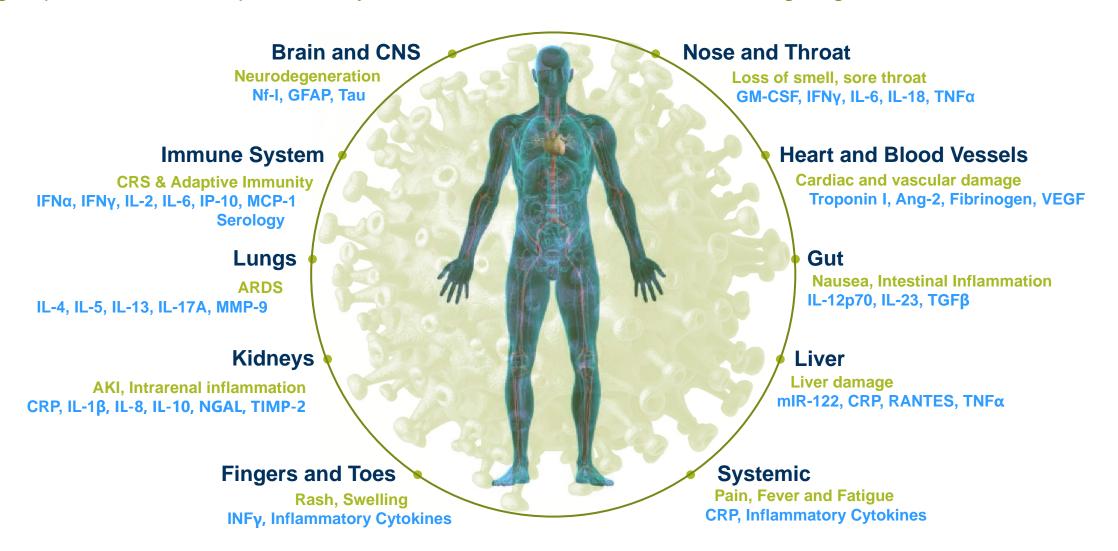
5

Comprehensive High-Definition Kinetic Picture – PPH Ecosystem



SARS-CoV-2 Infection Impacts Multiple Organ Systems

Digital protein biomarkers provide early detection of acute disease and monitoring longitudinal manifestations



Diagnostic Entry Accelerated by COVID & RADx

ACCELERATORS



Expanding Mfg and building Capacity, Regulatory & Quality Capabilities

Established **FDA** Credibility & Enabling major **Payor + Population Surveillance Studies**

Builds distributed infrastructure to attract and support future **Neuro CDx**

COVID PPH Think Tank connecting dots for Nextgen viruses and drug & vaccine efficacy

GO-TO-MARKET



Strategic focus on standing up at **University / Payors** with On-Site CLIA labs

Focus **Pre & Asymptomatic testing & Expand** to Nasal/Oral Swab, Saliva, DBS

Value Proposition - Accuracy, Price, High-Throughput, independent supply chain and Home Collection and Pooling

Liquid Biopsy expanding into Proteomic

Prostate Cancer



Only assay capable of precise measurement of PSA following prostatectomy





Al Roker



Pancreatic Cancer



LIF levels with Simoa showed elevation in Pancreatic cancer patients

Better correlation with response than FDA-approved CA19-9 marker

Alex Trebek



Brain Cancer



Neurofilament Light Chain as Biomarker for Brain Metastases



Bo Biden



Breast Cancer



Protein Biomarkers
Measurement in
CancerSEEK yielding
False Negative Results

Simoa's ~700x lower LLOQ holds the potential for accurate early detection

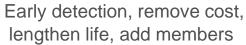
Angelina Jolie





Millions of Members







Inflammation

Infectious disease

Neurology

Cardio

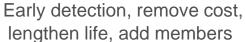






Millions of Members







Inflammation

Infectious disease

Neurology

Cardio

Massive Opportunity to Transform Healthcare

Five IRB's Underway



Known Biomarkers and Therapy

DEPLOY NOW

MS, Diabetes, COVID







Millions of Members



Early detection, remove cost, lengthen life, add members



Cardio

Neurology

Massive Opportunity to Transform Healthcare

Five IRB's Underway



Known Biomarkers and Therapy

MS, Diabetes, COVID

DEPLOY NOW





Known Biomarkers, No Therapy

FIND THERAPIES

Alzheimer's, Parkinson's, ALS











Millions of Members



Early detection, remove cost, lengthen life, add members



Massive Opportunity to Transform Healthcare

Five IRB's Underway



Known Biomarkers and Therapy

DEPLOY NOW

MS, Diabetes, COVID





Known Biomarkers,
No Therapy
FIND THERAPIES

Alzheimer's, Parkinson's, ALS









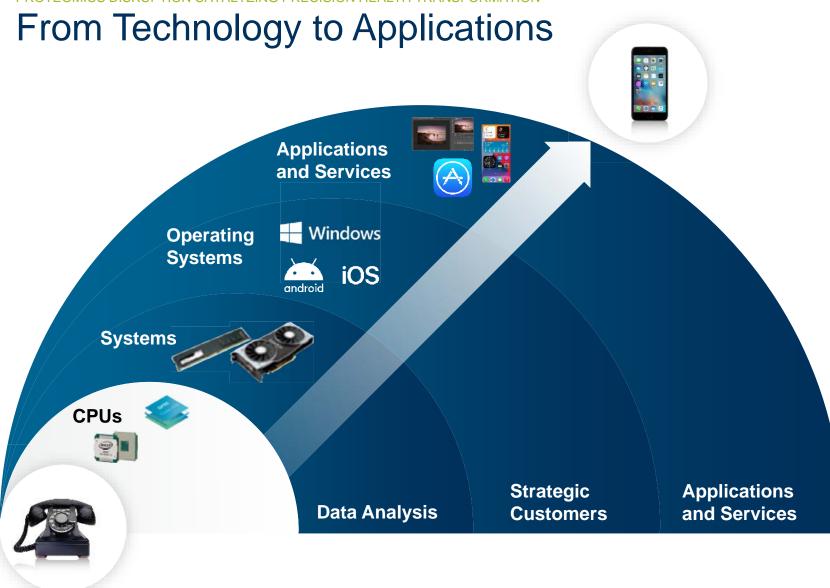


Known Therapy, No Biomarkers

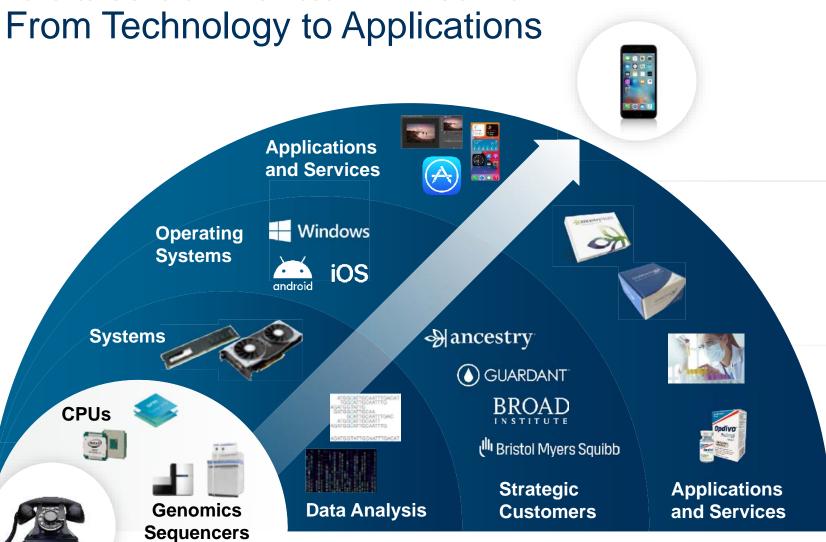
BIOMARKER DISCOVERY

Liquid Biopsies, Long-term Surveillance Study









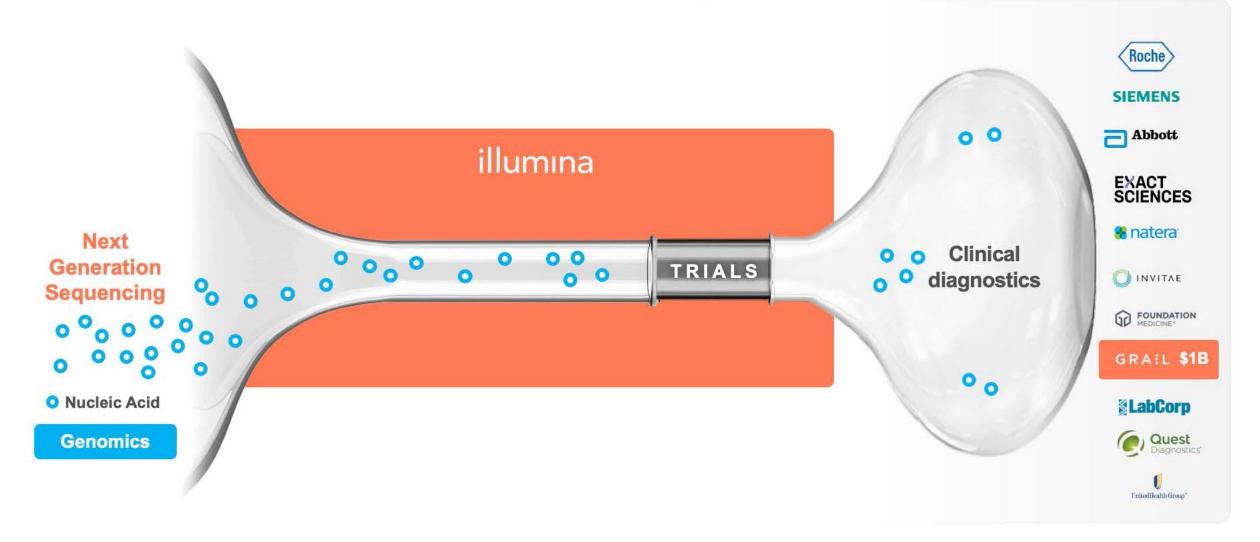
\$16B VALUE ACQUIRED Thrive \$2.5B

Illumina \$48B VALUE ACQUIRED Grail \$8B From Technology to Applications **Applications** and Services **Windows Operating Systems** iOS Systems **∌** ancestry (GUARDANT BROAD **CPUs** 9 Bristol Myers Squibb **Applications Strategic Genomics Data Analysis** and Services **Customers** Sequencers **Proteomics**

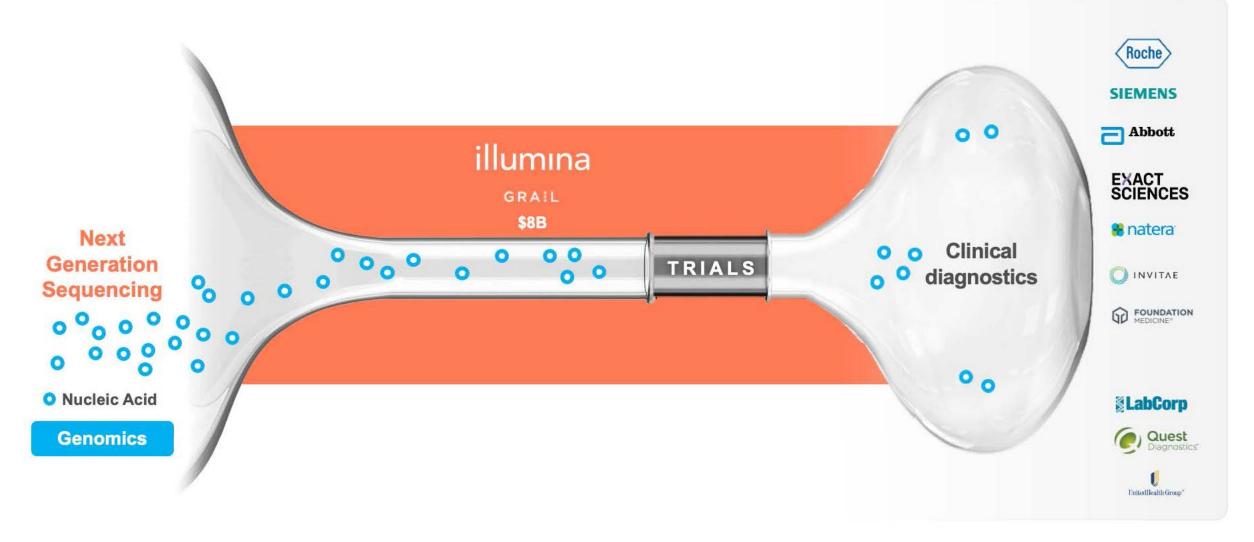
Exact Sciences \$16B VALUE **ACQUIRED** Thrive \$2.5B

> Illumina \$48B VALUE **ACQUIRED Grail \$8B**

Digital Molecules of Life; QTRX Focused on Proteins & Sensitivity



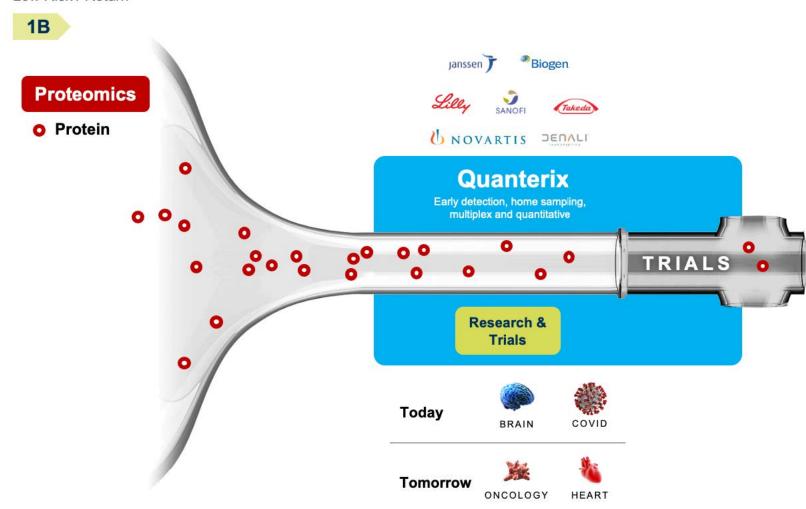
Digital Molecules of Life; QTRX Focused on Proteins & Sensitivity



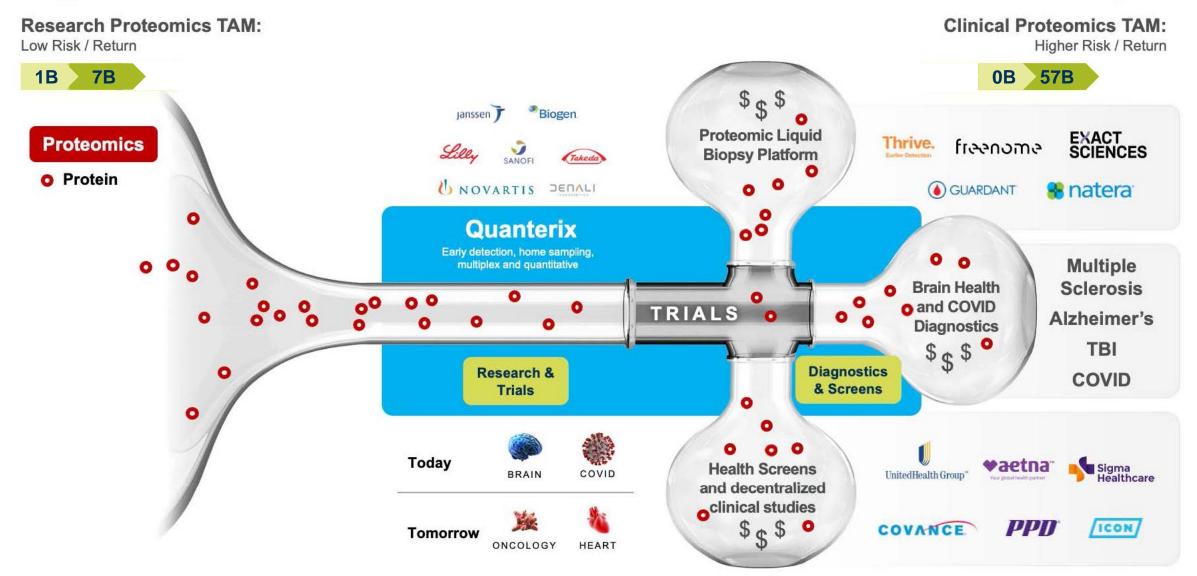
Digital Molecules of Life – QTRX Focused on Proteins & Sensitivity

Research Proteomics TAM:

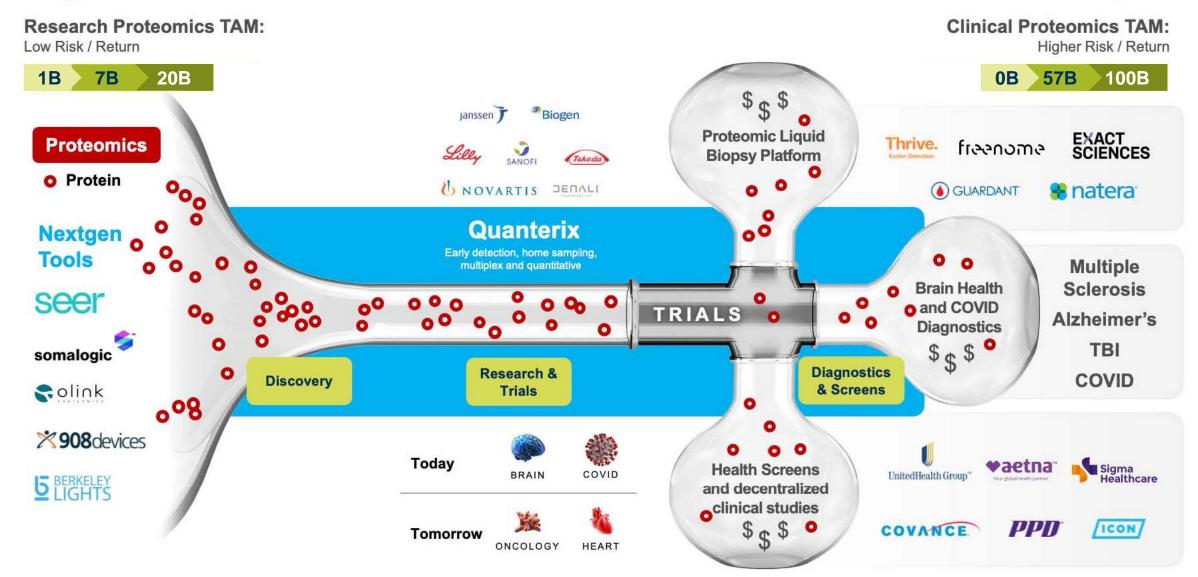
Low Risk / Return



Digital Molecules of Life – QTRX Focused on Proteins & Sensitivity



Digital Molecules of Life – QTRX Focused on Proteins & Sensitivity



Compelling Thesis to Lead Proteomics into Precision Health Era

Execution

2 to 3x Value Creation

RESEARCH & DISCOVERY

\$1B to \$20B TAM



COVID

RADx

- ✓ Validation: Pubs, Trials, Pharma Adoption
- 2014-20 Rev: \$0 \$85m; 55% CAGR* Public
- Raised \$350m; Valuation \$30m to \$2.0 B
- ST & LT Growth Catalysts
 - Increase penetration Neuro, COVID, Immuno
 - Menu expansion & drug / population trials
 - Scale consumables growth New COO
 - Expand into upstream Discovery tools

* 2017-2020 3Yr CAGR

Aspiration

10 to 15x Value Creation

DIAGNOSTICS & HEALTH SCREENS

\$100B TAM



- Abbott, Siemens, Pharma Advances
- Payor Acceptance for Cost / Outcome
- NIH, RADx & FDA Ecosystem expansion
- **⊘** ST & LT Growth Catalysts
 - COVID EUA's
 - Neuro Tool Entry Leverage COVID
 - Liquid Biopsy Entry (Partner or Vertical)
 - Advance Payor "Outcome Leapfrog"



Aspiration 2.0

DIGITAL HEALTH





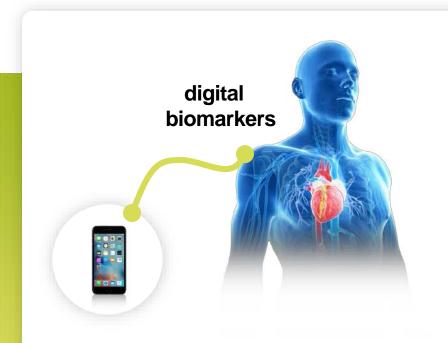














Biogen to Launch Pioneering Study to Develop Digital Biomarkers of Cognitive Health Using Apple Watch and iPhone

BIOMARKER WATCH Fitbit on Steroids!!



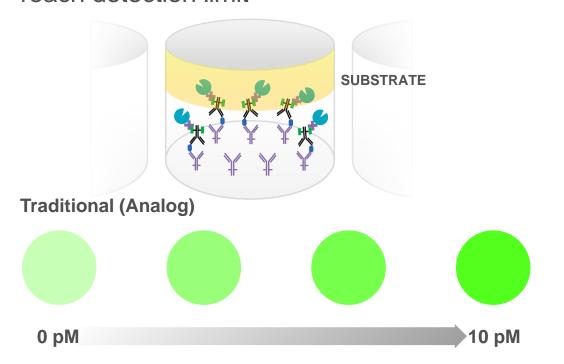
Appendix

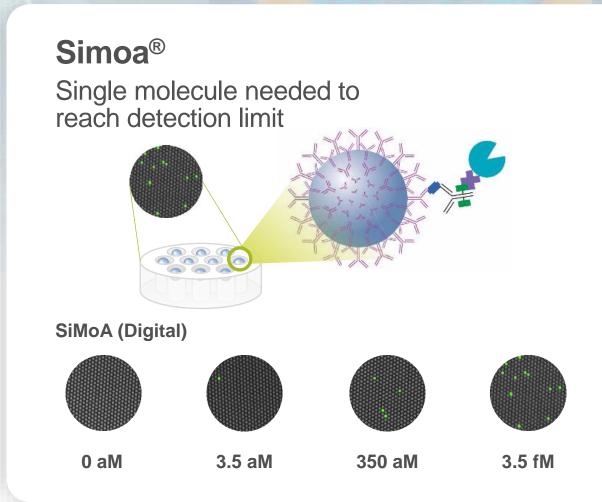


Our Single Molecule Arrays (Simoa®) technology allows for the ultimate sensitivity by detecting a single molecule

Traditional ELISA assay

Millions of molecules needed to reach detection limit





Quanterix Product Offerings

INSTRUMENTS

Bead based





List Price: \$200k 311 Installed (149 HD-X, 162 HD-1)

HD-X

Fully automated

Neuro Focused

SR-X

List Price: \$75k 175 Installed Benchtop Neuro Focused

Planar



SP-X

List Price: \$75k 35 Installed Benchtop Onco Focused

ASSAY KITS





Consumables & Reagents

400+ assays Homebrew kits Singleplex and Multiplex

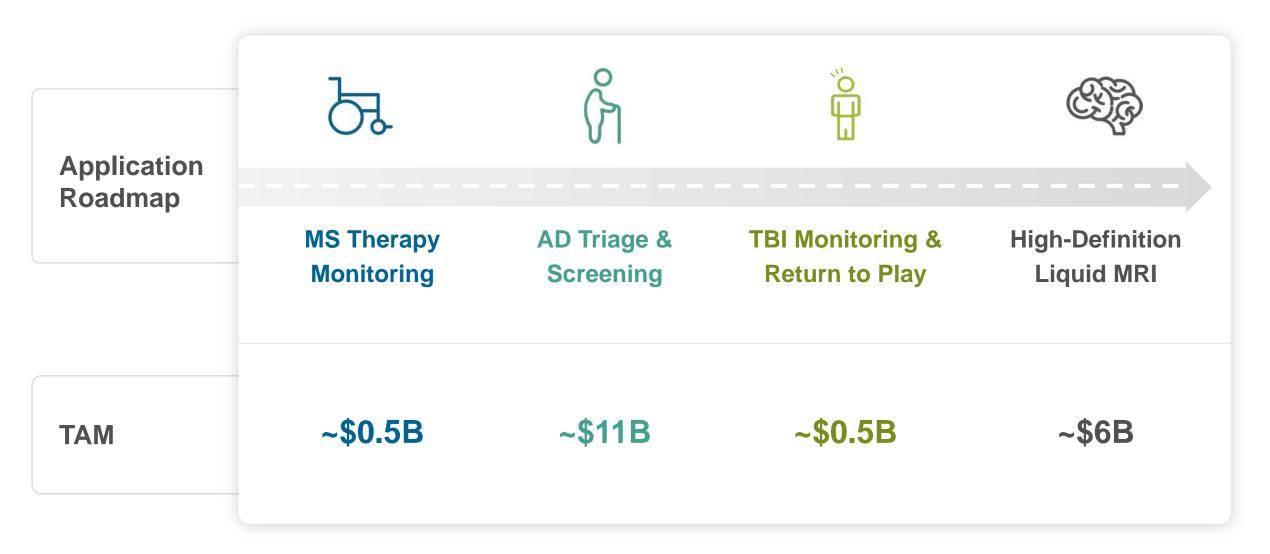
SERVICES



Accelerator

Contract research Custom assay development & reagent production CLIA and LDT capabilities Drug Rescue

Brain Diagnostic Health Roadmap



SIMOA supporting Leading-Edge COVID Research

Keys to developing Ultra-Sensitivity driven differentiated offering Long-Term COVID impact

Measuring changes in viral load in response to therapeutic mAb & convalescent plasma

Measuring relationship between adaptive immune response, viral load, and clinical outcome

Establishing biomarker profile (serology, cytokines, cardiac, neuro) in Long-haulers

Assessing impact of emerging COVID19 variants on accuracy of molecular and antigen tests

Longitudinal Surveillance testing for COVID19 population prevalence

Measuring Type I Interferon response in prediction of disease severity in COVID19

Measuring biomarkers of axonal injury and neurodegeneration in COVID patients

Early detection of cytokine dysregulation (alternatives to anti-IL-6 therapies in cytokine release)

Profiling serum biomarkers for associations between neuro. & pulmonary disease in Long-haulers

Characterization the role of autoantibodies against Type I IFN in severe COVID-19 patients



























There are major LB applications across the disease continuum

Neuro Diagnostics Market anticipated to also expand across these 5 Application Vectors . . .

<u>Applications</u>	Early detection*	Diagnostic aid and risk stratification	MRD* / active surveillance (AS)	Therapy guidance	Therapy monitoring
	Unaffected patients		Affected patients		
<u>Description</u> •	Enable early cancer detection in asymptomatic patients	 Complement existing tests to improve diagnosis or risk stratification 	 Measure MRD* and intervention response and longitudinally monitor for recurrence 	 Inform tx. selection in advanced disease for targeted therapy or IO 	 Monitor response or resistance to therapy in advanced disease
<u>Current state</u> <u>of market</u>	Nascent outside of CRC	Niche and nascent	Deployed in clinical trials	Becoming established clinically	Nascent beyond EGFR T790M in NSCLC
Addressable patients (U.S.)	Tens of millions (Eligible adults)	Millions (High risk adults)	Millions (Early stage patients)	Hundreds of thousands (Late stage patients)	
Market leaders / front runners	EXACT SCIENCES	✓ veracyte.	natera* Conceive. Deliver. Thrive.	GUARDANT	Roche
	GRA!L	MD*Health. biotechne	GUARDANT	FOUNDATION MEDICINE	QIAGEN

Disruptive Opportunities with High-Growth potential



Opportunity

Value of ultra-sensitivity

Initial Quanterix partnerships

TAM

Population Health

Assessment of large, population-level datasets

- Ability to assess previously undetectable biomarkers
- Unlock visibility into earlier disease states
- Non-invasive sample compatibility (low volume)

Major US Payor group

~\$5B



Decentralized (virtual) Trials

At-home assessment of subject health status

- Non-invasive and self-sampling compatibility (e.g., finger prick, dried blood spots)
- More frequent & robust Longitudinal Data
- Higher Subject Compliance & throughput (Enhanced Speed & Lower Cost)
- Inclusion of Under-Served Populations

Major CRO Labs

> **RADx Funding**

~\$5B

