

J.P. Morgan Healthcare Conference 2022

Kevin Hrusovsky, Chairman & CEO January 13, 2022

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. Forward-looking statements in this presentation are based on Quanterix' expectations and assumptions as of the date of this presentation. 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 presentation 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 forward-looking statements contained herein to reflect any change in expectations, even as new information becomes available.



Transforming Medicine with Hyper Sensitive & Quantitative Low Plex NextGen Proteomics



Precision Health Vision

2021 Achievements



Symbiotic Relationship

COVID - Neuro & RUO - Diagnostics



Alzheimer's Opportunity

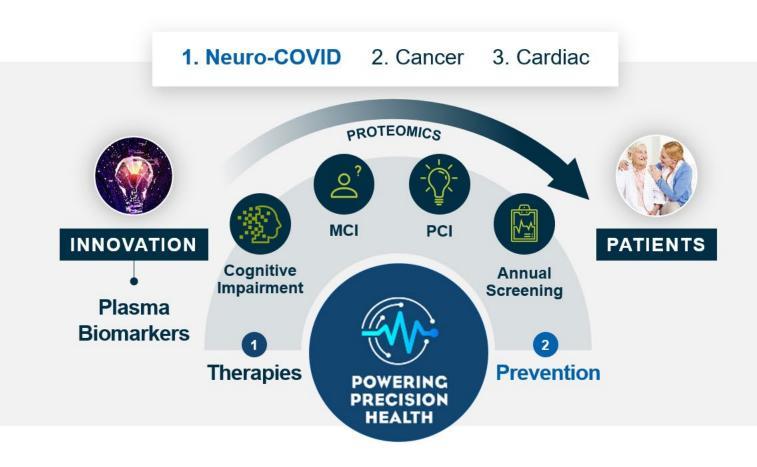
2022 Objectives

Investment Thesis

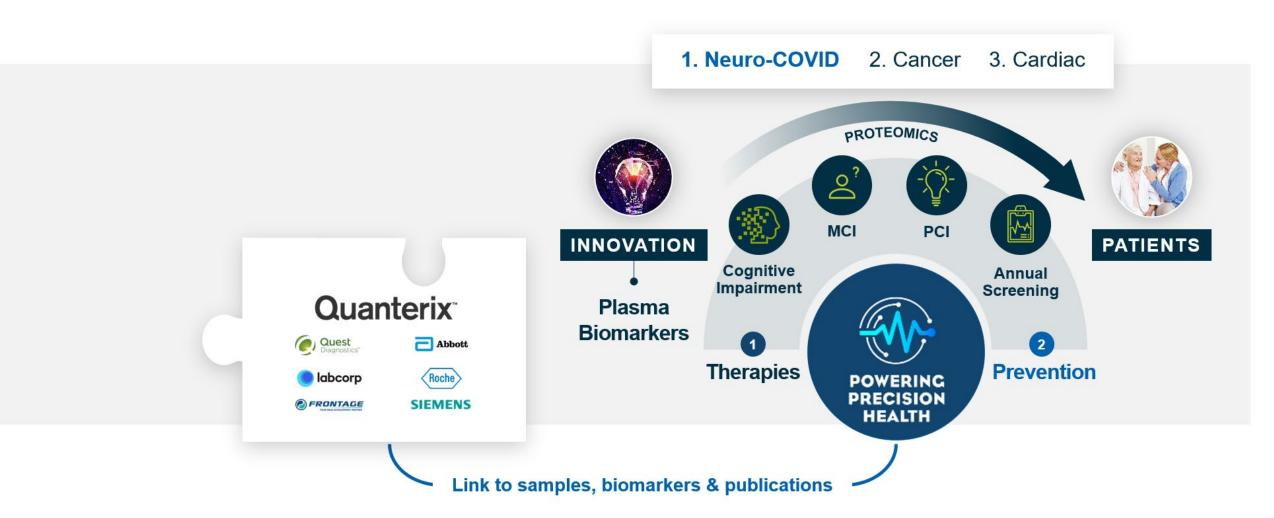


Questions & Discussion

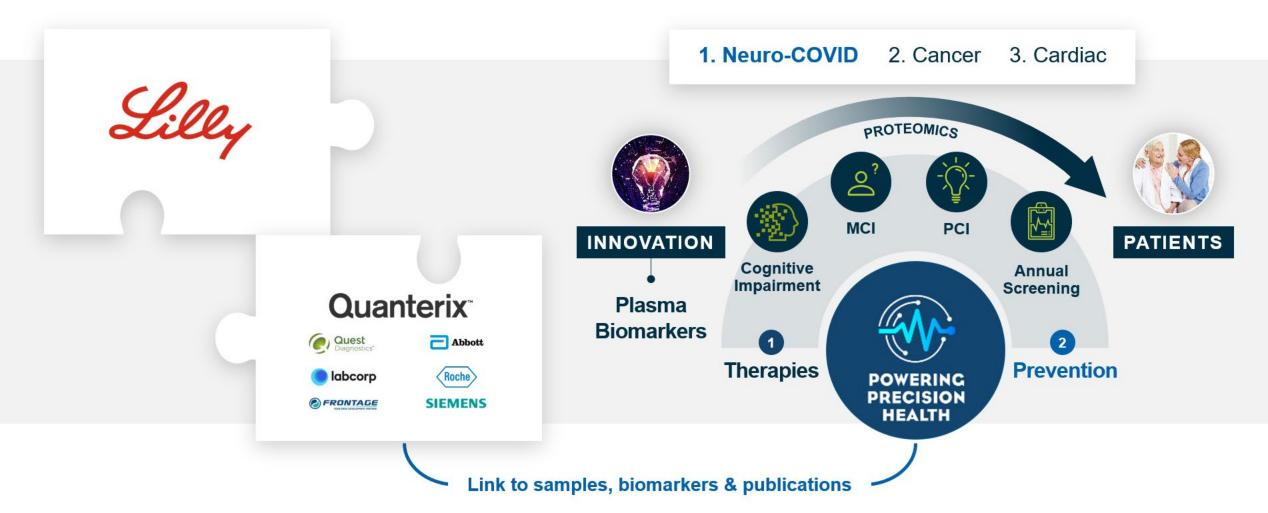




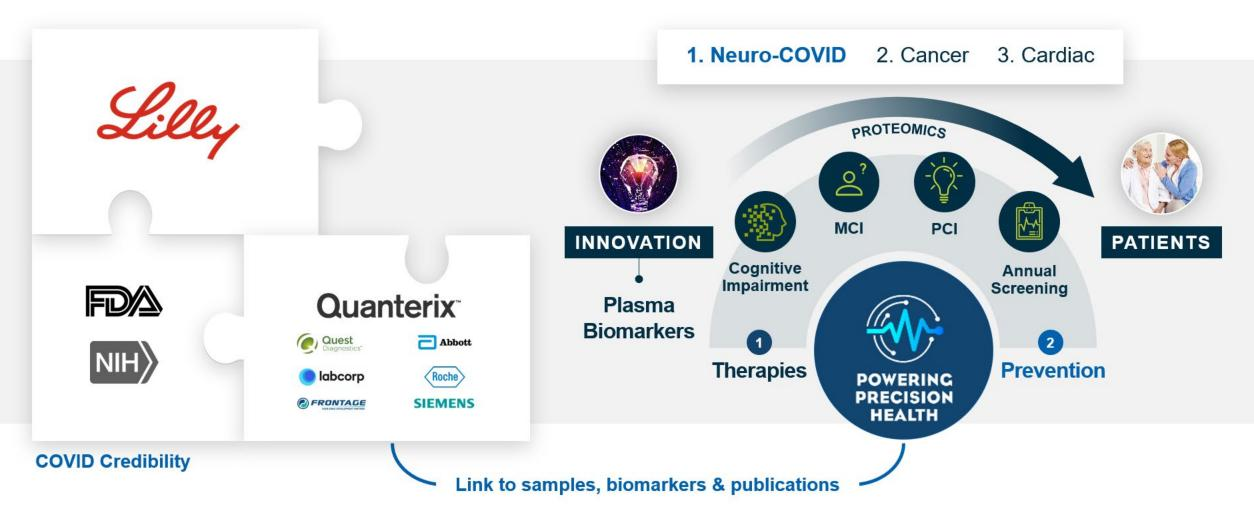




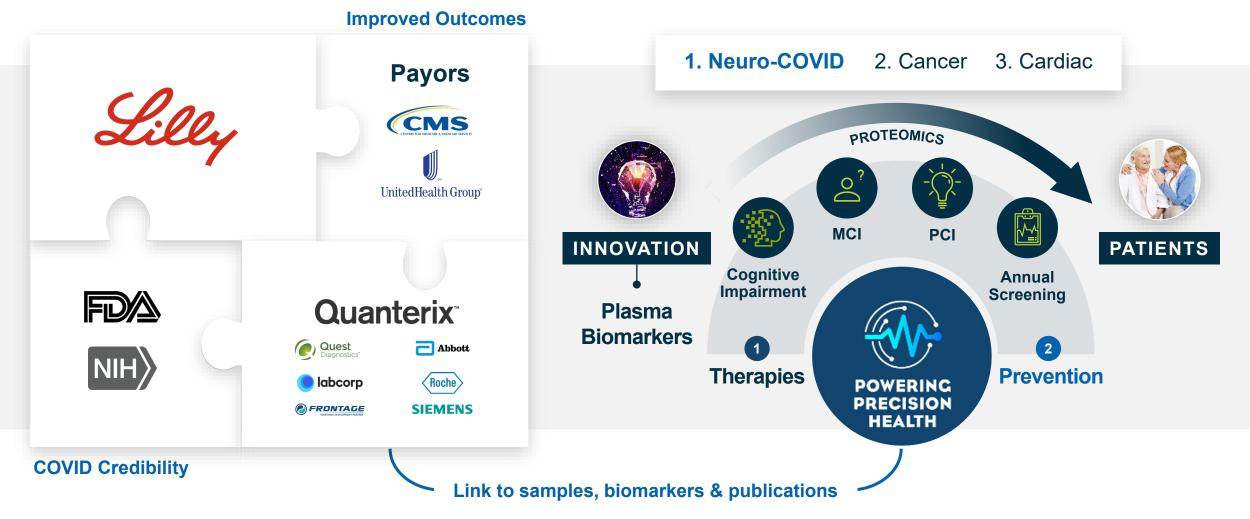














Blood-based biomarkers

are fast becoming gold standard surrogates for traditional Tau or Abeta PET imaging

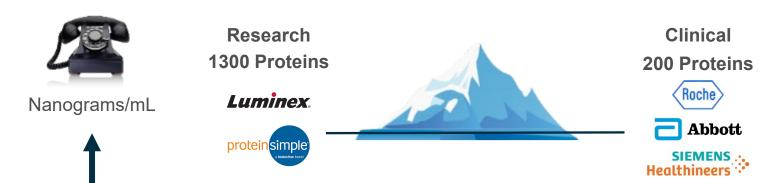


Ever since the FDA approved Biogen and Eisai's Alzheimer's disease treatment Aduhelm, the buzzword among competing companies has been "biomarker." And that's just what Eli Lilly has to offer with two new analyses on their candidate donanemab.

> - Fierce Biotech, "Lilly adds on to biomarker bombardment in Alzheimer's with 2 new donanemab analyses"



Digital Biomarkers Sensitivity Unlocking Proteomics



LATE: INVASIVE

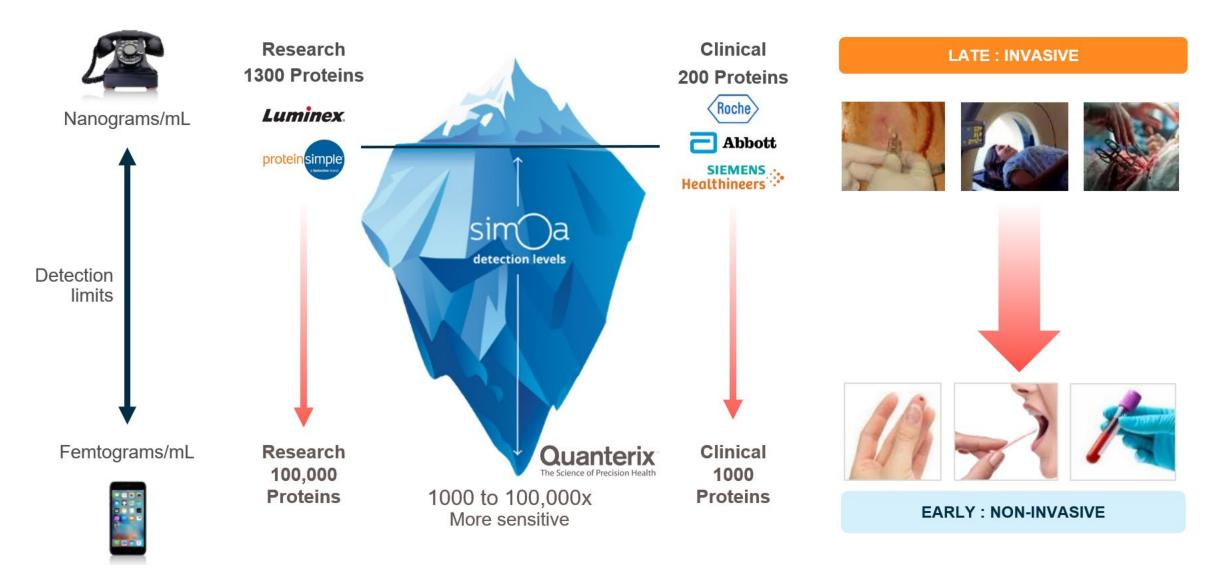




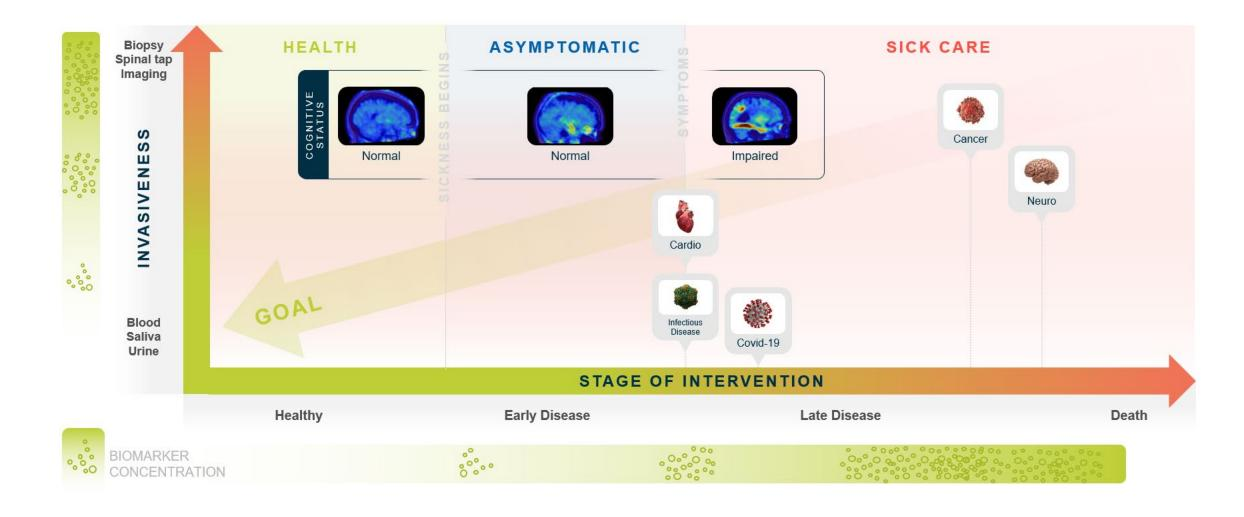




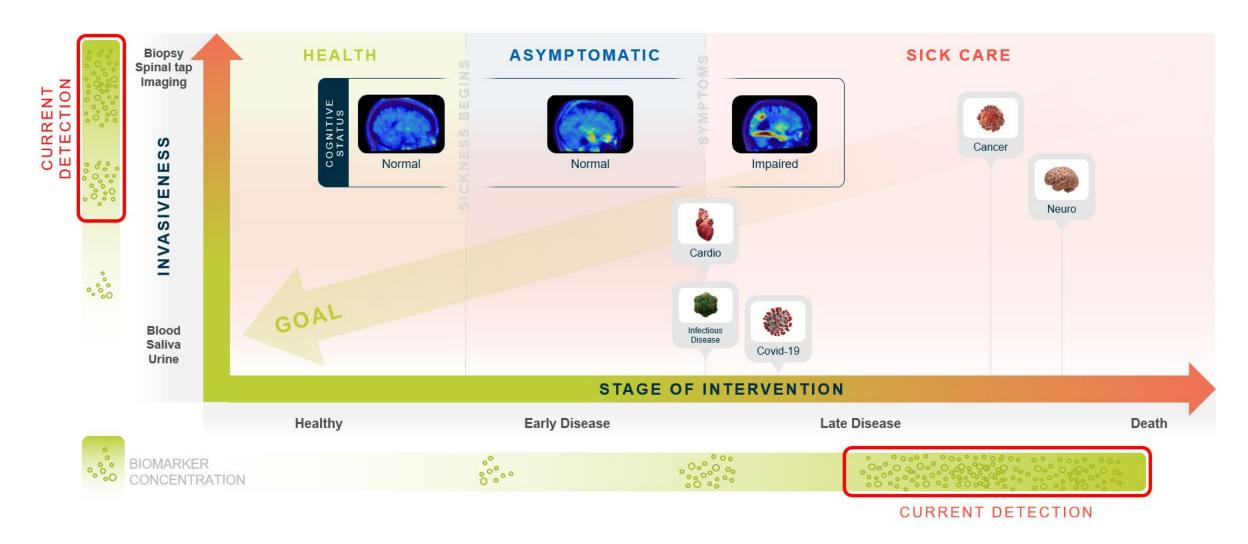
Digital Biomarkers Sensitivity Unlocking Proteomics



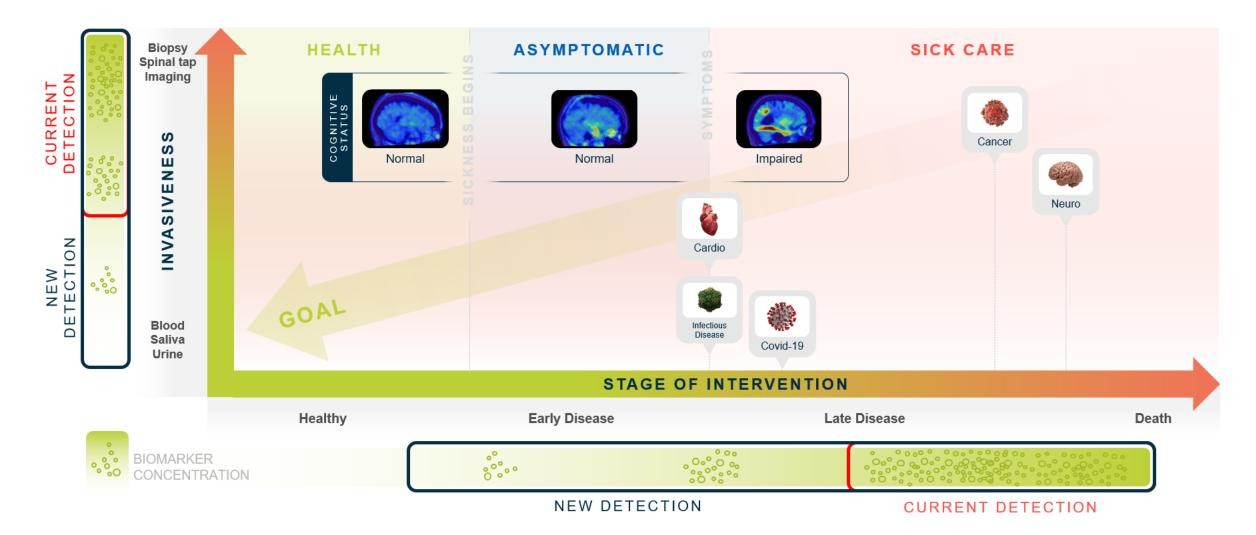




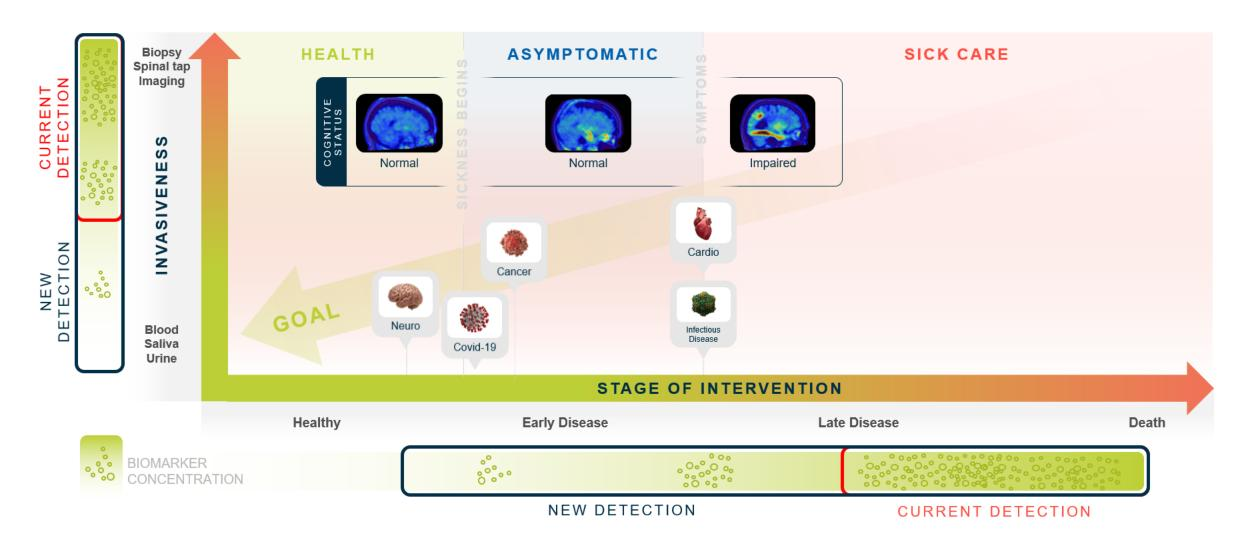














Proteomics Value Chain

Shift From Symptomatic To Asymptomatic

	Discovery	TRANS Research	LATIONAL Drug Trials	Dx & Screens
Plexity	100+	1-50	1-10	1-10
TAM 2025+	\$6 Bn	\$2 Bn	\$2Bn to \$12 Bn Increased adoption	\$30Bn to \$100 Bn Asymptomatic + Payors
Cost & Thru-put	Cost			Thru-put Low



Proteomics Value Chain

+ Relevant Proteins

Shift From Symptomatic To Asymptomatic **TRANSLATIONAL Discovery** Research **Drug Trials Dx & Screens Plexity** 100+ 1-50 1-10 1-10 \$2Bn to \$12 Bn \$30Bn to \$100 Bn TAM \$6 Bn \$2 Bn 2025+ **Increased adoption Asymptomatic + Payors** High Cost & Thru-put Thru-put Cost Low NextGen Tools **Quanterix Today** Abbott **SIEMENS** Market Quanterix 2025+ + Early Detection + Home, less invasive + Early Detection

+ Early Detection

+ + Quantitative / Precise

+ Multiplex

+ Quantitative / Precise

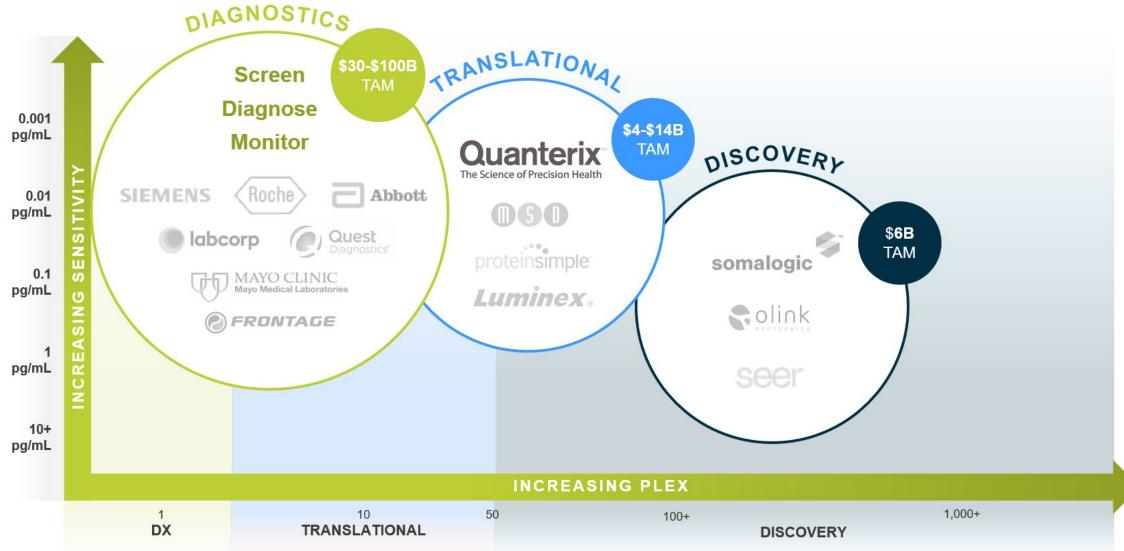


Sensitivity

+ Quantitative / Precise

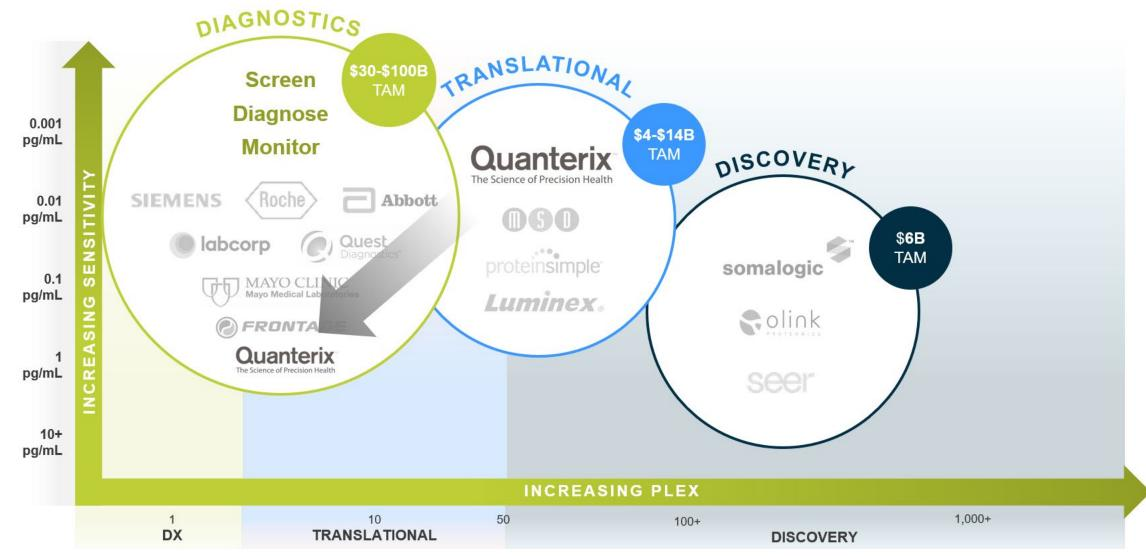
+ Home, Less Invasive

Proteomics Landscape – Sensitivity – Multiplex play a key role

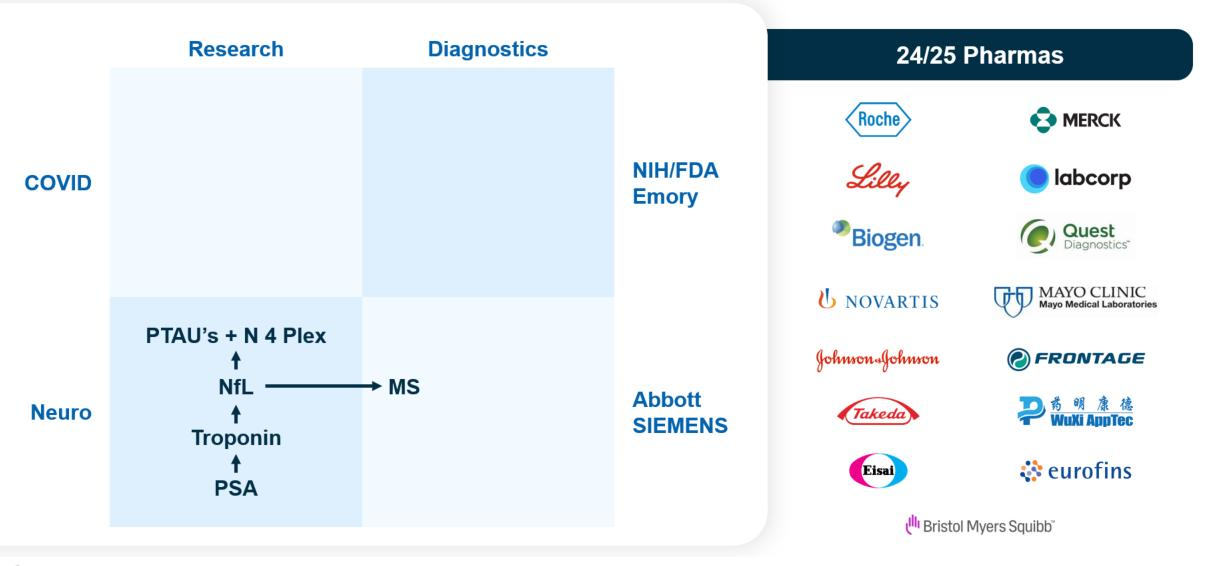




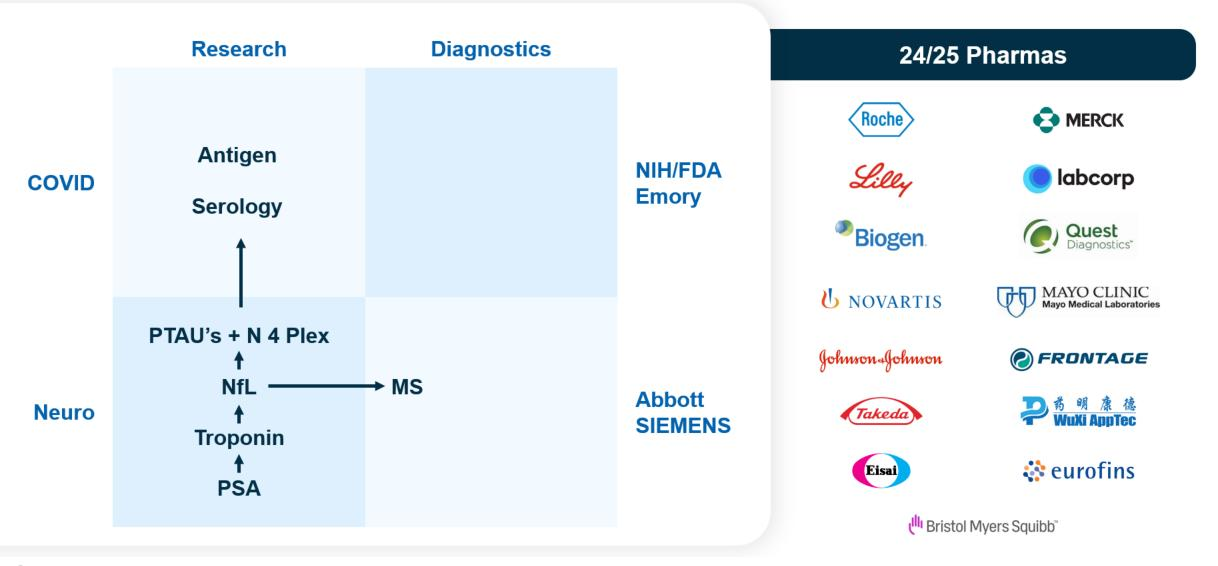
Proteomics Landscape – Sensitivity – Multiplex play a key role



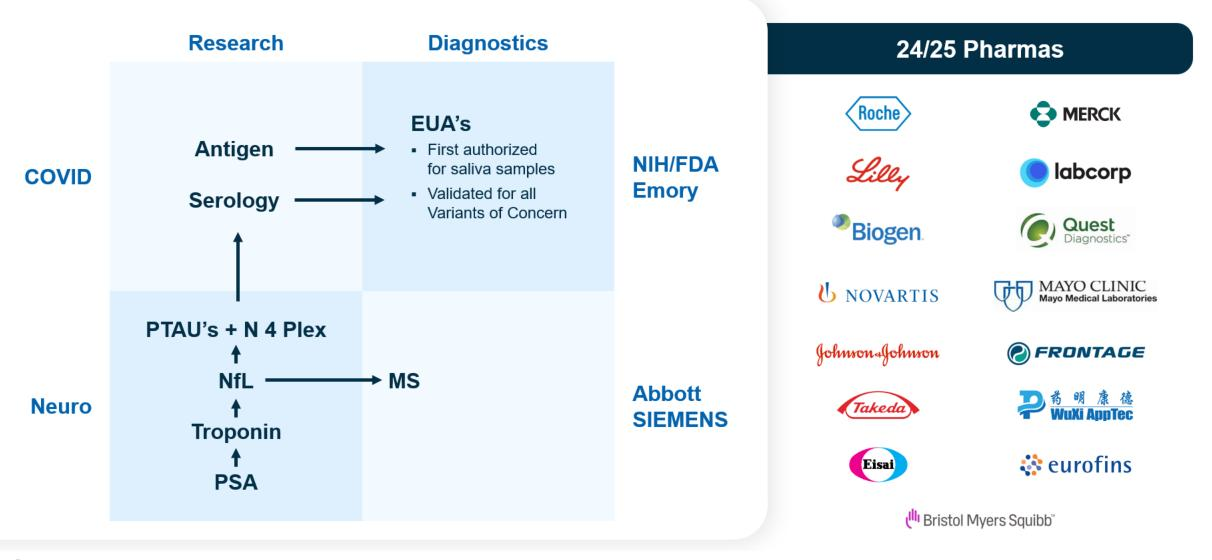




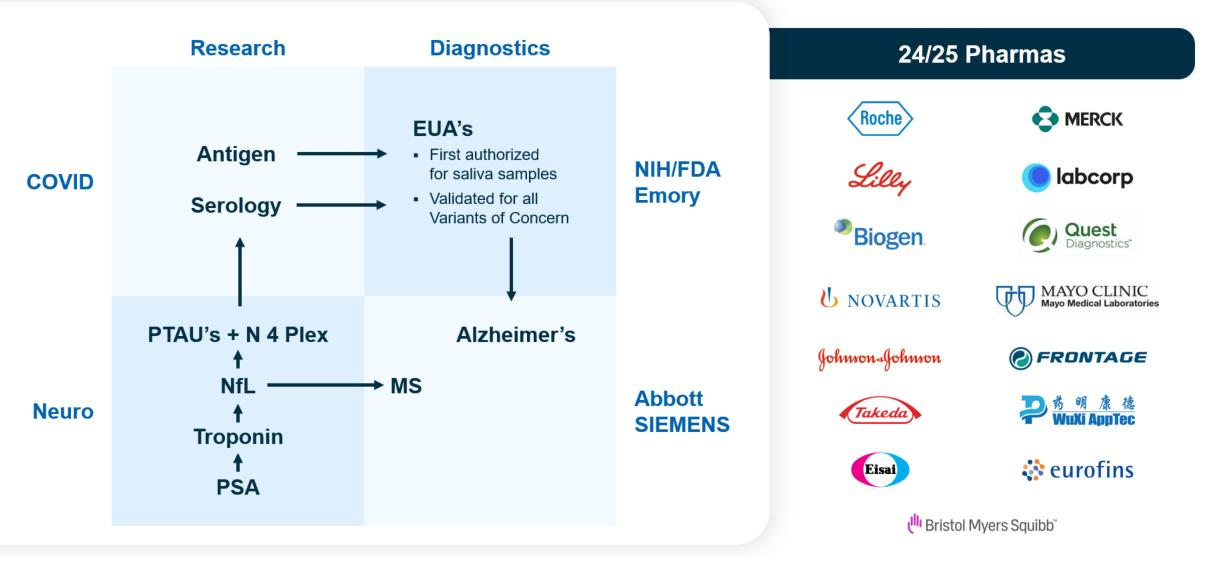




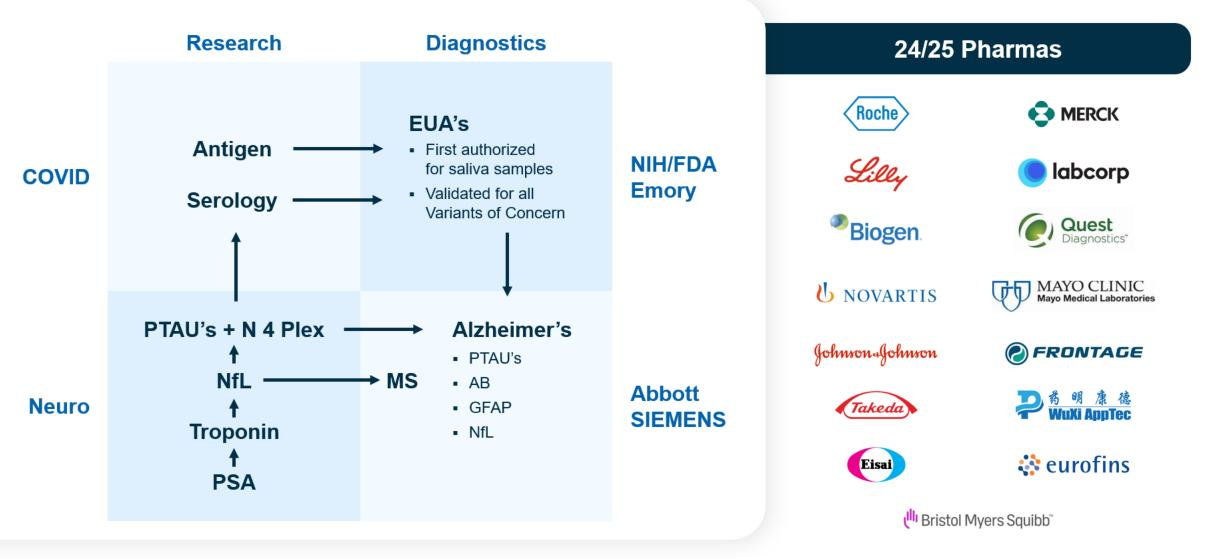




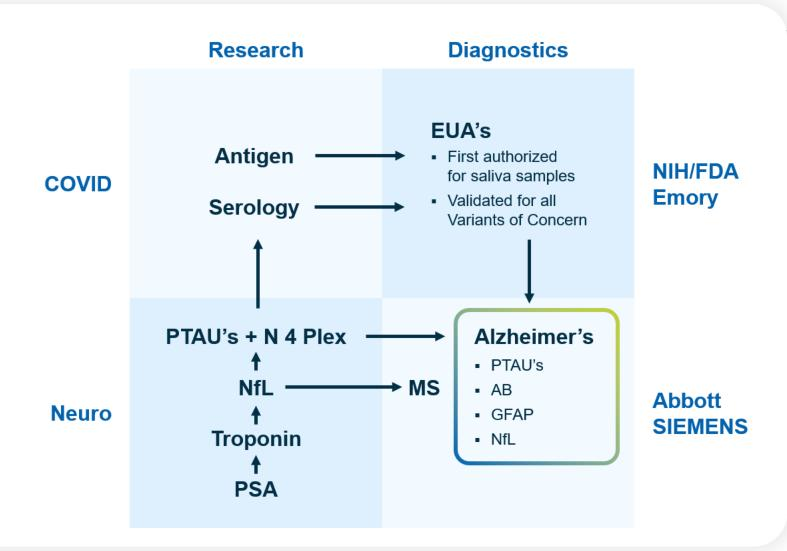












24/25 Pharmas





























Bristol Myers Squibb والله



Our 2021 Results

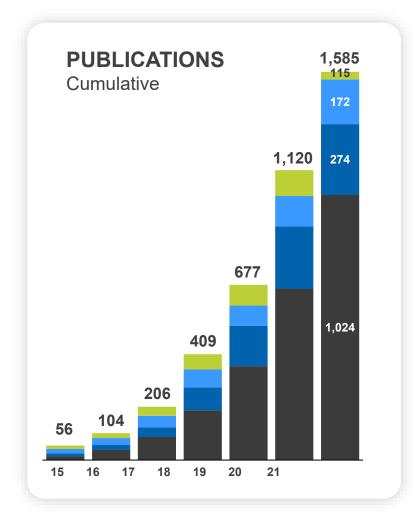
Continue superb execution & value creation via commercializing disruptive innovation

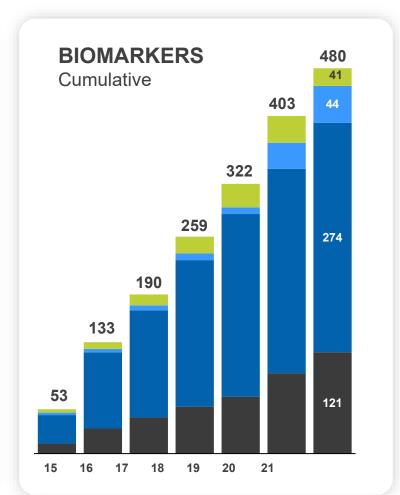
FY 2021 — RESULTS		
30-40% CAGR in RUO Product (consumables & instruments) revenue increase of 84% yoy	②	
Achieved plasma pTau 181 FDA Breakthrough Designation for Alzheimer's diagnostic / triaging	②	
Record 465 new publications Academic pull through brings SIMOA publications to 1,585	Ø	
Instrument install base grows 32% to 708; exceeding HDX target of 105 placements by 28 Shift to HDX vs. SRX provides up to 3x pull through		
Lilly and Biogen employ SIMOA for blood plasma biomarker detection in Trials Rapidly expanding RUO collaborations for DX entry	Ø	
Strengthened balance sheet raising 287M, \$410M on balance sheet end of Q3. RUO crossover 4Q'23	Ø	
100x Sensitivity Demonstrated in the Field	Ø	

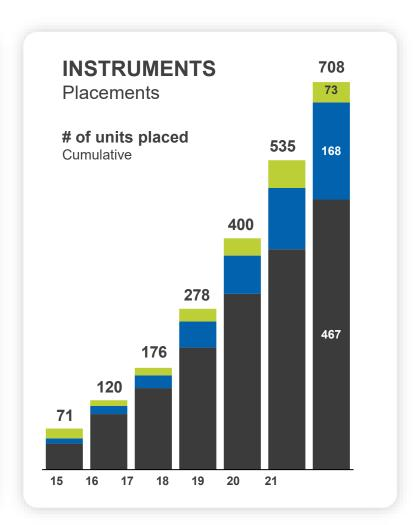


Scientific Validation Driving Adoption

2021 Advances







Infectious Diseases

Others

Immunology & Oncology

Neurology

^{*} Q4 2021 and FY 2021 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, 2021 are ongoing and could result in changes to such estimated information.



Accelerators Driving Rapid Adoption for Transforming Drug Development

Research Institutions



Biopharma



CROs & Other





Superb Execution Continues

Strategic Growth Indicators

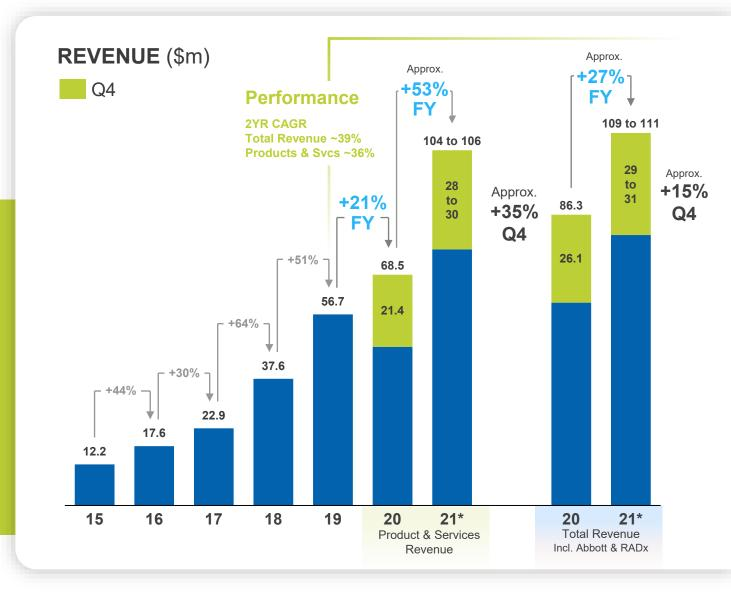
Approximate*

100% Record Annual Consumables revenue growth, driven by strong utilization

Proportion of HD-Fleet on new 68% HD-X technology at YE2021

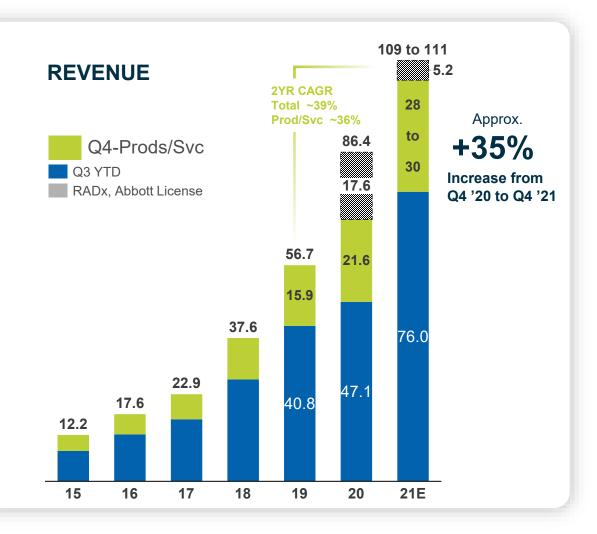
Instruments installed, with 133 215 HD-Xs (74% New) and 78 SR-Xs & SP-Xs

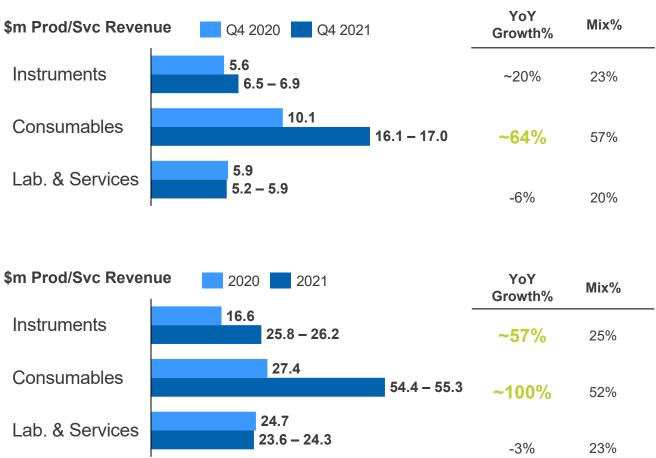
^{*} Q4 2021 and FY 2021 information are preliminary, unaudited estimates and are subject to the Company's normal quarter and year-end close procedures. These procedures and the audit of the Company's financial statements for the year ended December 31, 2021 are ongoing and could result in changes to such estimated information.





Q4 & FY2021 – Revenue Growth







Compelling Thesis to Lead Proteomics into Precision Health Era

Execution

2 to 3x Value Creation

TRANSLATIONAL

\$4B to \$14B TAM²



- ∀alidation pTau181 Breakthrough Device Designation
- ✓ Strategy: Fortify Moat w/ 100x sensitivity for Neuro
- **⊘** Growth Catalysts
 - pTau's & Neuro Plex: AD Diagnostic Therapies
 - Academic publications
 - Revitalized Neuro drug trials
 - Scale, HDx Clinical, Accelerator Expansion & 100x

* 2017-2020 3Yr CAGR

. 409A Valuation YE 2014 2.

Based on independent third-party research report 2021



Compelling Thesis to Lead Proteomics into Precision Health Era

Execution

2 to 3x Value Creation

TRANSLATIONAL

\$4B to \$14B TAM²



- ✓ Validation pTau181 Breakthrough Device Designation
- Strategy: Fortify Moat w/ 100x sensitivity for Neuro
- **⊘** Growth Catalysts
 - pTau's & Neuro Plex: AD Diagnostic Therapies
 - Academic publications
 - Revitalized Neuro drug trials
 - Scale, HDx Clinical, Accelerator Expansion & 100x

Aspiration

10 to 15x Value Creation

Dx & SCREENS

\$30B to \$100B TAM²



- Mounting Evidence − AAIC plasma pTau's
- Strategy: Single site IVD for NfL-MS / pTau-AD
- Longer-term Growth Catalysts
 - AD triage, screens, diagnostics & monitoring
 - Increase Accelerator LDT footprint and capability
 - Payor leapfrog for health screens

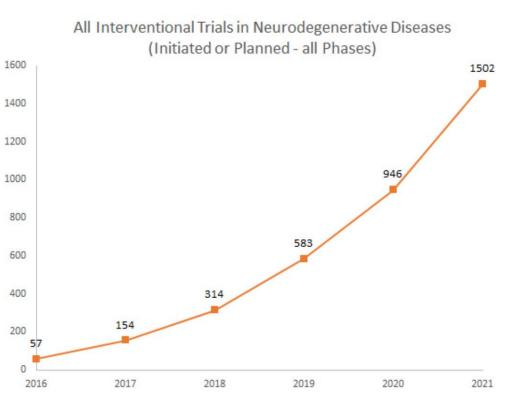
1. 409A Valuation YE 2014

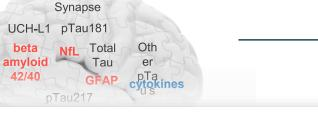
Based on independent third-party research report 2021

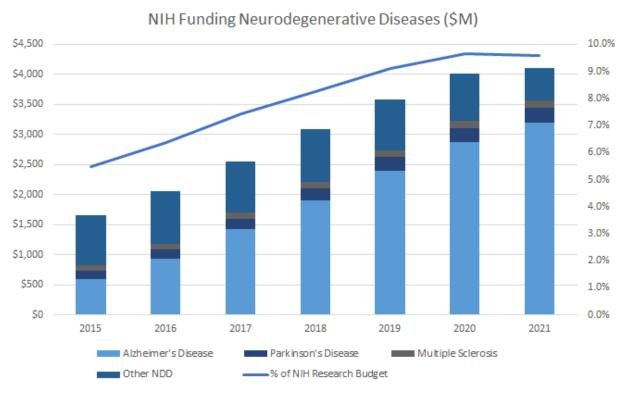


^{* 2017-2020 3}Yr CAGR

Entering the Neuro Decade..... with a Leadership Position







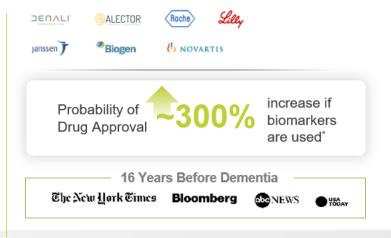


Alzheimer's poised for Transformation

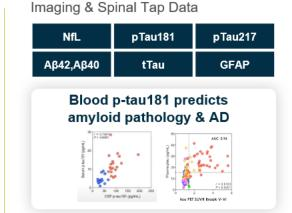
Drug-trials center-piece of near-term focus







2 Clinical Utility



Clinical Trials
FDA Breakthrough Designation
10% market share = \$50M

Source: Bio Industry Analysis; Clinical Development Success Rates (June 2016) LINK



Alzheimer's poised for Transformation

Drug-trials center-piece of near-term focus

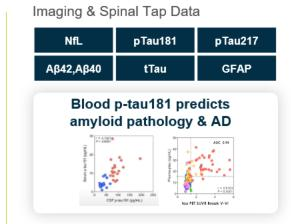




TAM <\$0.5B



Clinical Utility



3 Dx & Screens







- Multiplex panels >90% AUC
- 5.7M AD in US to grow to 13M by 2050
- Early screening key to deliver real therapeutic effect
- Global cost of AD is over \$1T

Clinical Trials

FDA Breakthrough Designation

10% market share = \$50M

Enter Dx

FDA single site IVD approval

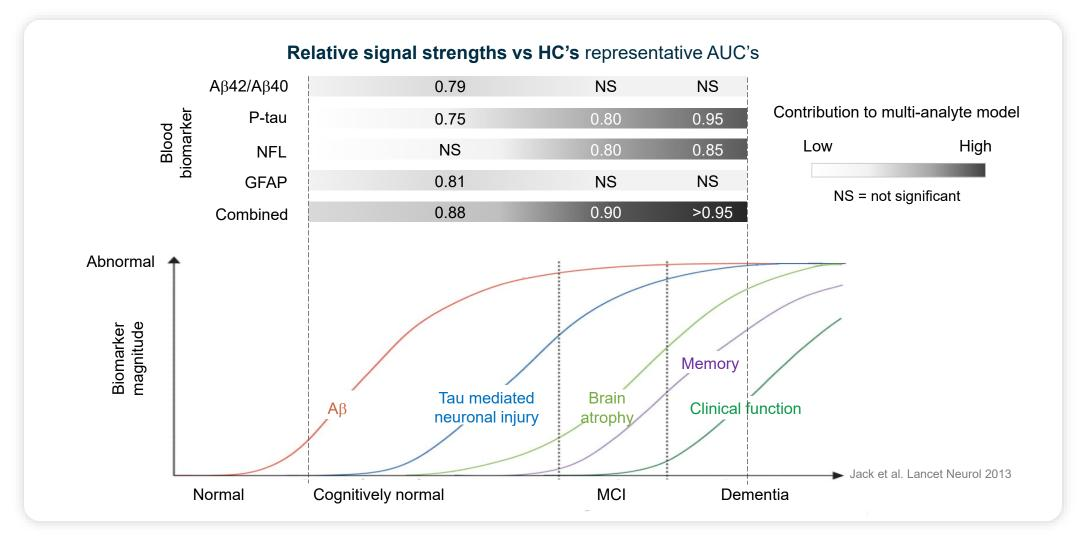
10% of this market share \$1B

Source: Bio Industry Analysis; Clinical Development Success Rates (June 2016) LINK



Putting it all together: multi-analyte approach to early detection

Multiple analytes are more accurate than any single analyte





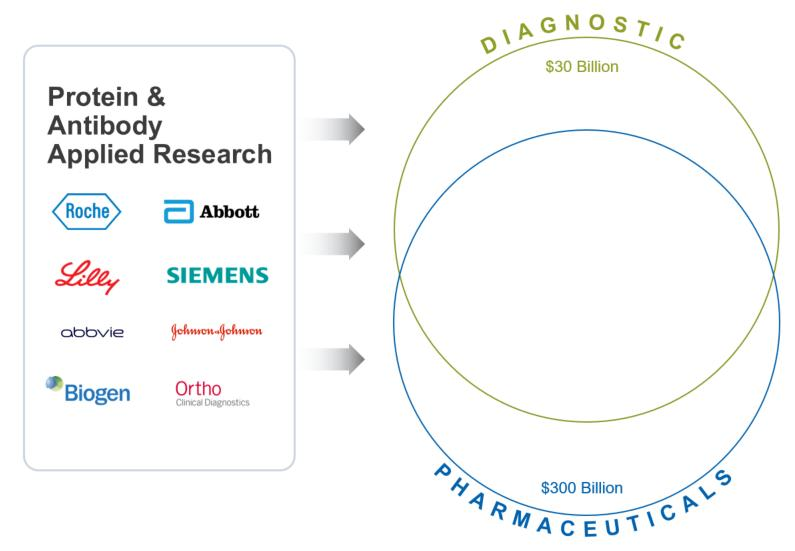
Blood Neurodegeneration Biomarkers: Quanterix is Clear Leader

Quanterix leads all other new proteomics platform entrants in its inclusion in published, peer-reviewed scientific literature and studies

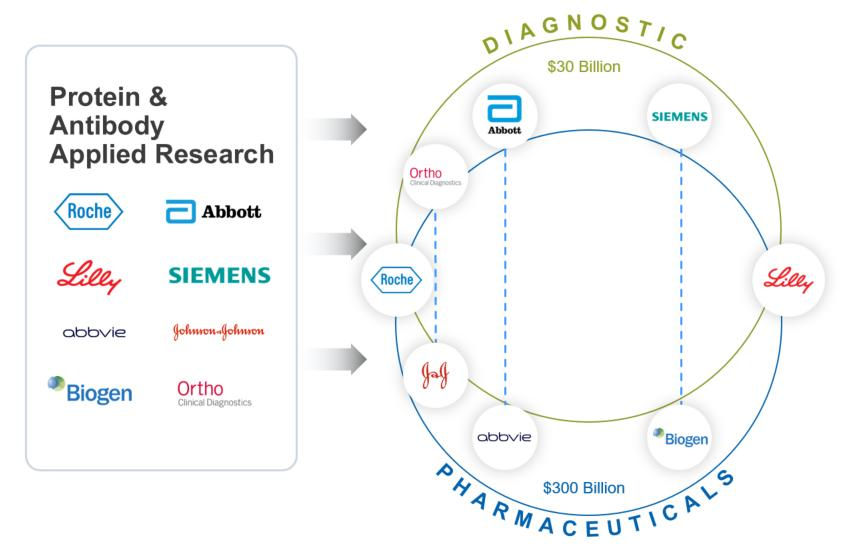
Alzheimer's	Parkinson's	ALS	MS	Concussion
	Alzheimer's Alzheimer's	Alzheimer's Parkinson's Alzheimer's Parkinson's Alzheimer's Parkinson's	Alzheimer's Parkinson's ALS ALS ALS ALS ALS	Alzheimer's Parkinson's ALS MS Alsheimer's ALS MS Alsheimer's MS Alsheimer'

1585
Publications

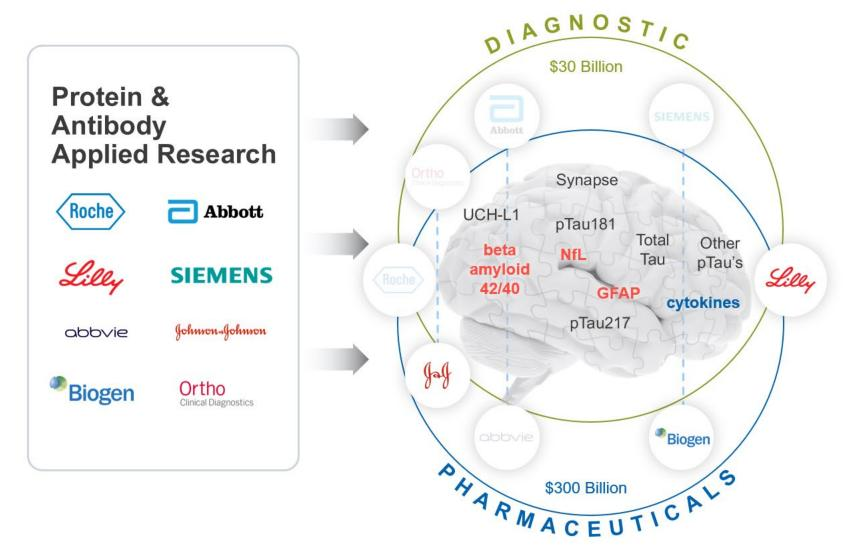




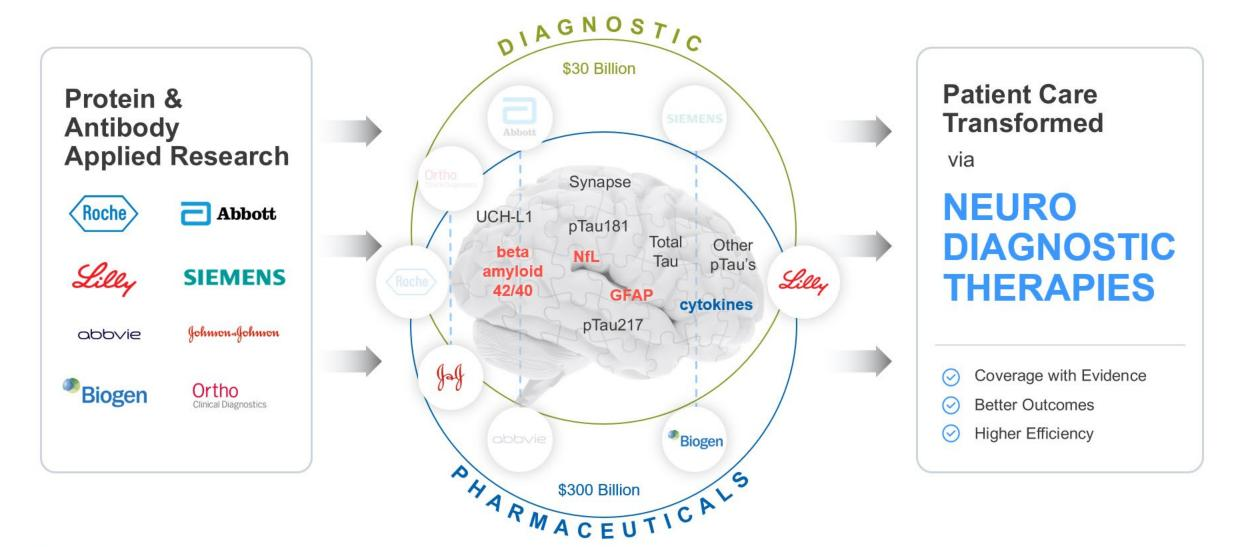










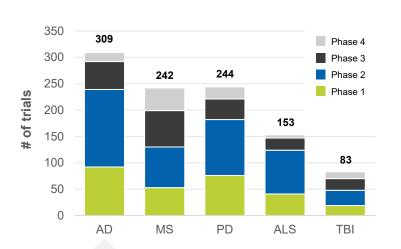


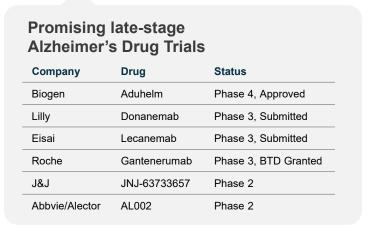


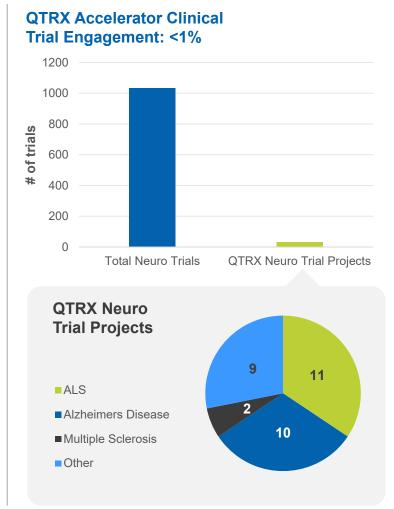
QTRX Neuro Clinical Trial Engagement

Significant Opportunity to Expand Neuro Clinical Trial TAM Capture Through Accelerator & CROs

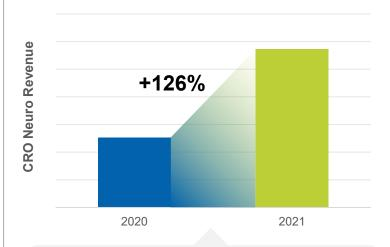
>1000 Active Clinical Trials in 2021







QTRX CRO Neuro Revenue (YTD Q3)

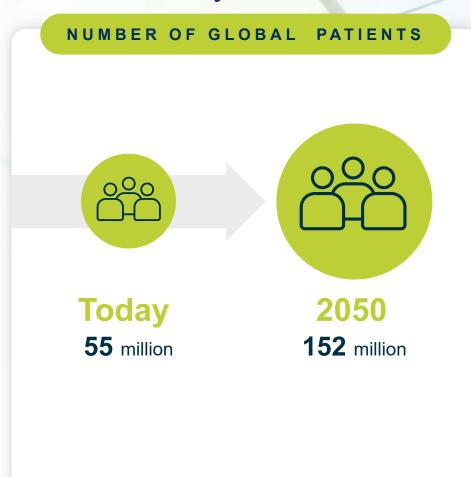


Highlights

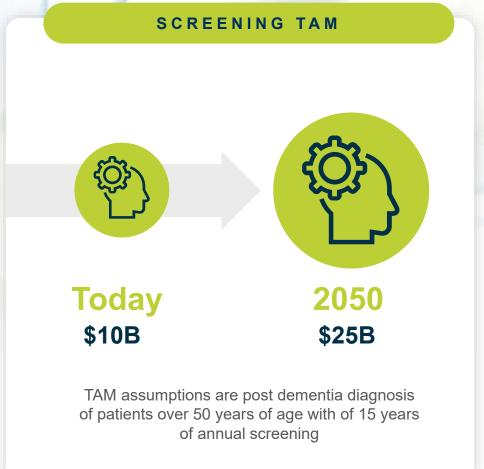
- 1000 active trials for AD, MS, PD, ALS, and TBI;
 QTRX participating in <1% through Accelerator
- Significant Opportunity to Expand Neuro Clinical Trial TAM Capture Through Accelerator & CROs



Worldwide Dementia Cases Expected to Triple by 2050 Assumes only 10% of the world has Diagnostics access





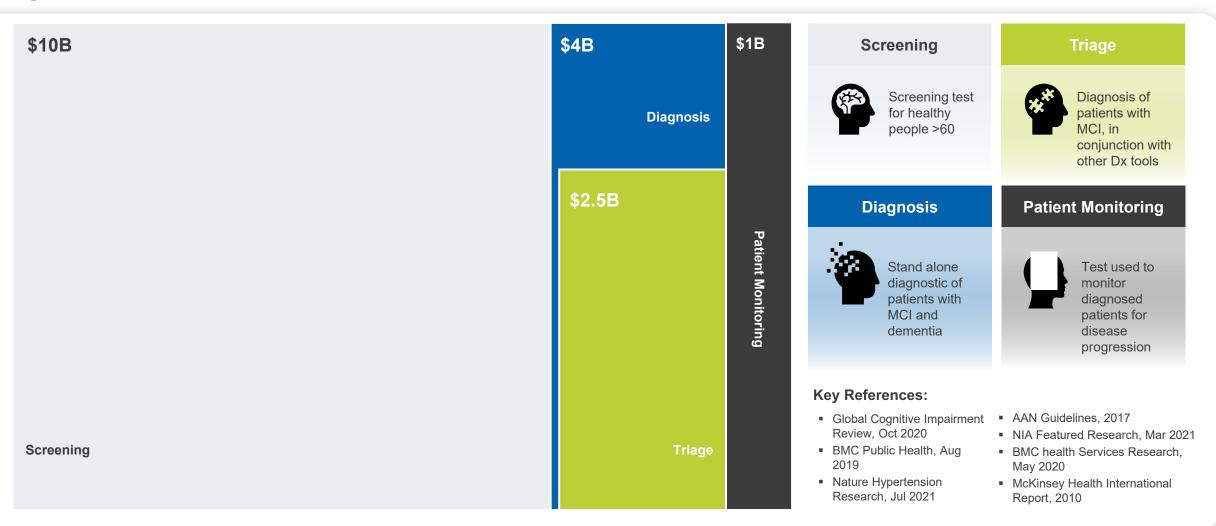


Source: www.alz.org



Alzheimer's Disease Diagnostic Testing Market: \$15B

Global





CMS Draft National Coverage Determination Guidance

Highlights the vital role for biomarker measurements in pre- and post-market clinical trials











Pharma Coverage with Evidence Development Enabled by Simoa Biomarker Testing:

Accelerated and economical Phase 4 trial enrollment through plasmabased biomarker screening

Quantitative drug efficacy evidence development via correlation with PET imaging & cognitive assessment in postmarket clinical trials Non-CNS contributions to amyloid in the blood compartment limit the AD vs healthy control differential to just 10-15% for blood-based amyloid assays¹, making pTau singleplex and multiplex assays a better choice for enabling clinical discrimination





Simoa biomarker testing enables Pharma partners to establish clinical evidence



¹Schindler et al., Neurology, 2019

Accelerating Clinical Market Entry Alzheimer

Pharma Partnerships has the potential to Accelerate and De-risk Market Penetration

Rule Out

Primary care setting

Identifies the 'worried well'

Maximize sensitivity



Maximize specificity

Neuro specialist

Identifies Alzheimer's patients

Clinical trials support both 'rule in' and 'rule out' use cases — each served by setting different clinical cut-offs

TRIAGE INITIALLY, SCREEN ULTIMATELY

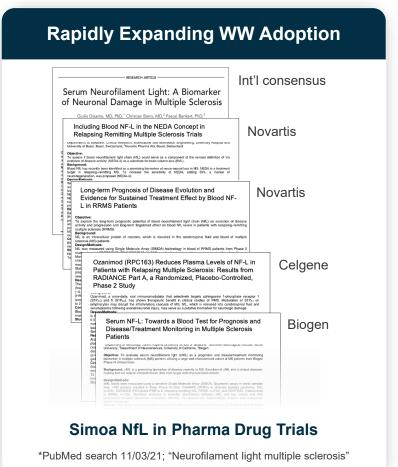
Stages	Timeline	Investment	TAM	
Laboratory Develop Test (multiplex / singleplex)	2023	\$10M-\$15M	Today	
Laboratory IVD (multiplex / singleplex)	2024	\$40M-\$60M	- \$15B NA	\$60B ROW
Distributed IVD (multiplex / singleplex)	2025	\$100M	2050 \$45B NA	\$180B ROW



Simoa NfL in Multiple Sclerosis

Std of Care for MS Monitoring \$3,000 MRI Simoa NFL blood test "MRI shows the ashes, NfL shows the match." David Leppert, M.D., Univ of Basel



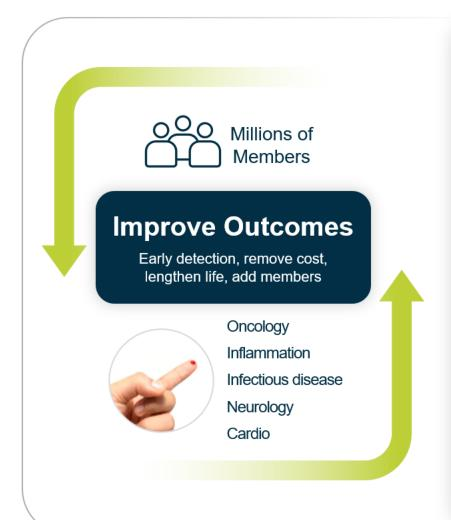


Source: PubMed









Opportunity to Transform Healthcare



Known Biomarkers, & Therapy

DEPLOY SHORT TERM

MS, Diabetes, COVID







Opportunity to Transform Healthcare



Known Biomarkers, & Therapy

DEPLOY SHORT TERM

MS, Diabetes, COVID





Known Biomarkers, No Therapy

FIND THERAPIES

Alzheimer's, Parkinson's, ALS













Opportunity to Transform Healthcare



Known Biomarkers, & Therapy

DEPLOY SHORT TERM

MS, Diabetes, COVID





Known Biomarkers, No Therapy

FIND THERAPIES

Alzheimer's, Parkinson's, ALS











Known Therapy, No Biomarkers

DISCOVERY

Liquid Biopsies, Long-term Surveillance Study



Our Objectives in 2022

Continue superb execution & value creation via commercializing disruptive innovation

FY 2022 — OBJECTIVES	
30-40% CAGR in RUO. Maintain or expand utilization	②
Neuro LDT validation for pTau181 and / or Nf-L for MS	
Start AD clinical trial for pTau181 & other AD markers (panel)	②
Scale supply and RUO growth as DX entry increases RUO differentiation; Consider new pTau's/panels	
Expand pharma partnerships with greater Neuro drug trial penetration, i.e., target 10% penetration in 2025	Ø
Strategy: Expand RUO plexity capability to 20 plex by '24 and bolster LDT / SSIVD penetration	②
Launch 100X Beta Program via Accelerator by YE'22	



Aspiration 2.0

DIGITAL HEALTH





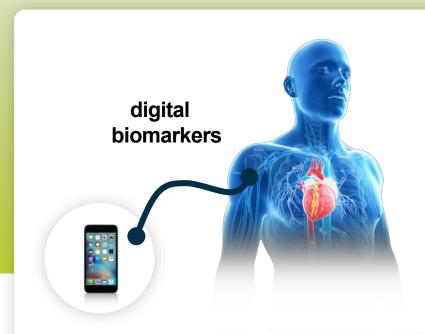












BIOMARKER WATCH Fitbit on Steroids!!





Poised to Disrupt Healthcare and Create Significant Value

EXECUTION: 2 - 3X VALUE CREATION

1 Unrivaled Sensitivity / Ecosystem

MARKET

- 2 Methodical market penetration strategy
 - 3 Proteomics; Better linked to Disease / Health

Quanterix
The Science of Precision Health

4 Validation: 19/20 top pharma, PPH, 1000+ trials

ASPIRATIONAL: 10 - 15X VALUE CREATION





Poised to Disrupt Healthcare and Create Significant Value

EXECUTION: 2 - 3X VALUE CREATION

1 Unrivaled Sensitivity / Ecosystem

MARKET

- 2 Methodical market penetration strategy
- 3 Proteomics; Better linked to Disease / Health

Quanterix
The Science of Precision Health

- 4 Validation: 19/20 top pharma, PPH, 1000+ trials
- Growth & Value; Razor razor blade, \$150M invested

PENETRATION

- 6 Low Risk / Solid Return + Uber Return Prospect
- 7 Track Record for Commercializing Disruption

ASPIRATIONAL: 10 - 15X VALUE CREATION

DIFFERENTIATOR

VALUE

Best in Class

Disrupt

\$1B to \$120B

New Answers



Holy Grail

1500+ pubs All Areas

Proven

Product Launches

Rapid Growth



Retail



Lynchpin

