

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 are based on Quanterix' expectations and assumptions as of the date hereof. 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 forward-looking 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



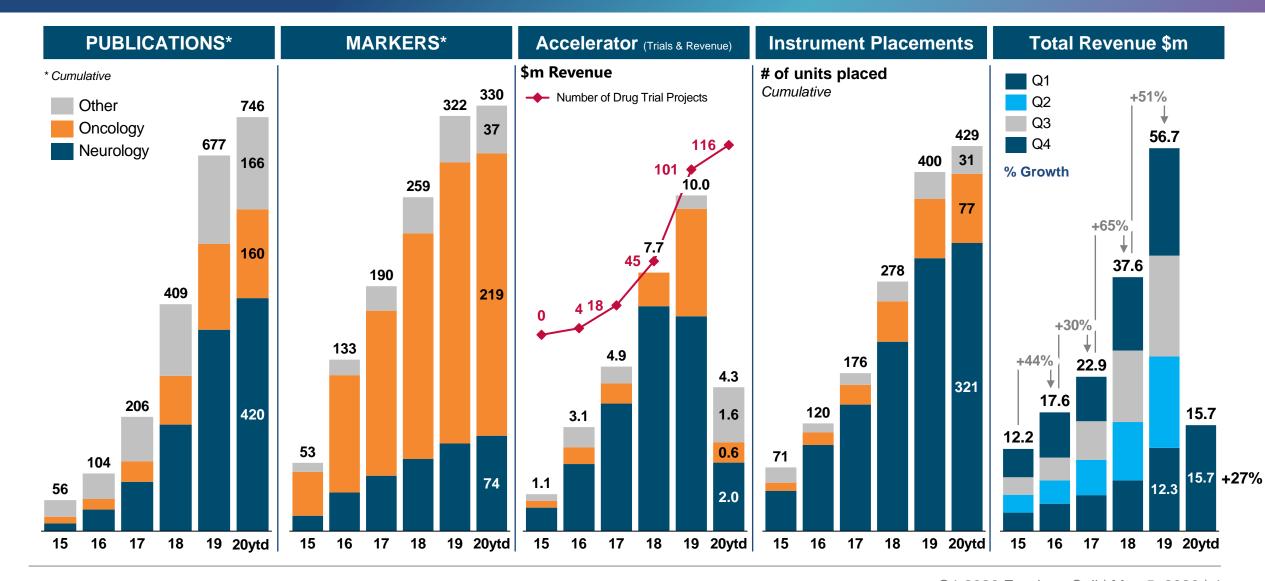
- I. Strategic and Financial Progress Kevin Hrusovsky Chairman, CEO
 - i. Q1 2020 Highlights
 - ii. Vision and Strategy
 - Executional / Aspirational value creation
 - Navigating COVID-19: Neuro + Near-term CAC PIVOT
 - iii. Beyond COVID-19

II. Financial Report – Amol Chaubal – CFO

III. Q&A

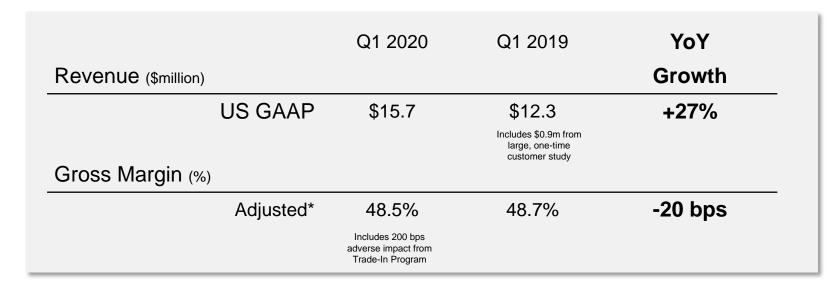
Continuing our journey into 2020, despite COVID-19 Headwind Further pivoting our focus to COVID/ID to drive breakout performance

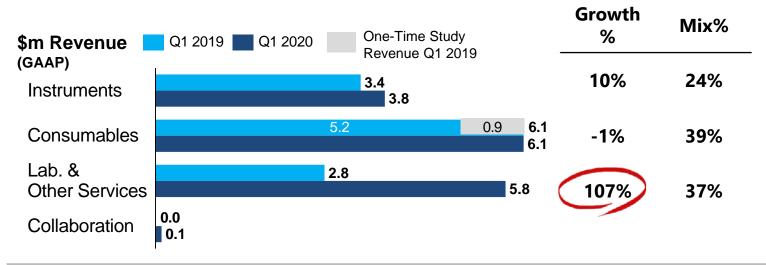




Q1 2020 Growth Led by Lab & Other services





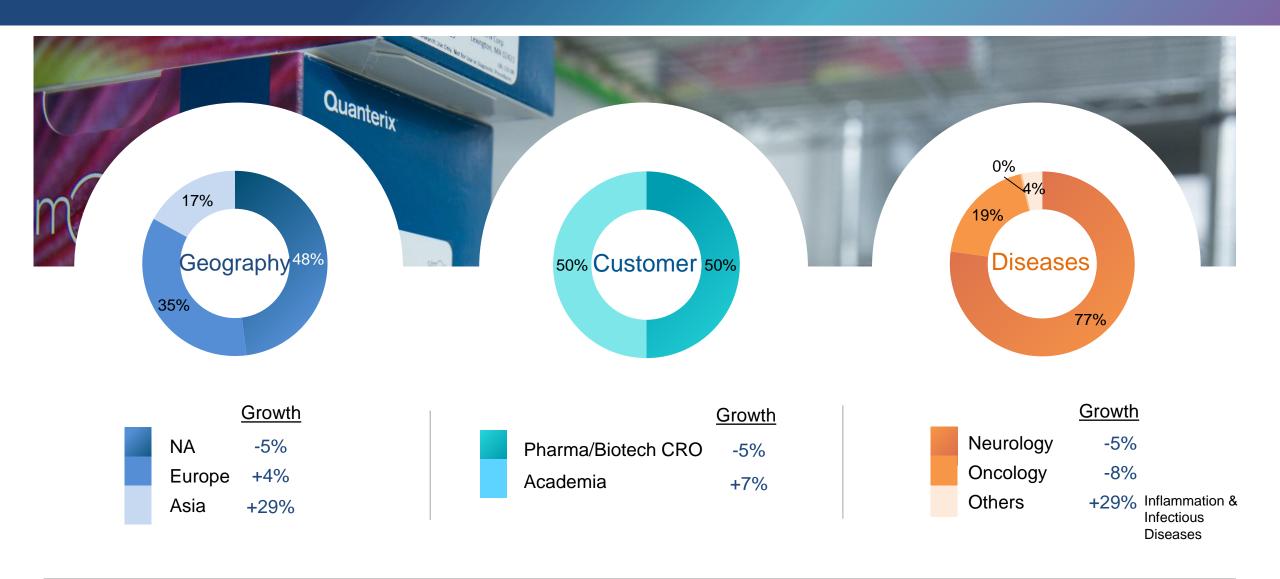


- Q1 Revenue growth +27%, despite install-base overhaul and COVID-19 impact
- Early steps taken to protect employees and implement resilience plan
- Accelerator capacity expanded to support customer disruptions & sustain clinical trials
- Engaged PPH collaborators worldwide to combat against COVID-19 pandemic
- Continued publication flow underscoring the growing relevance of our technology
- Reinforced the growing utility of Simoa in neurology research

Q1 2020 Earnings Call | May 5, 2020 | 5

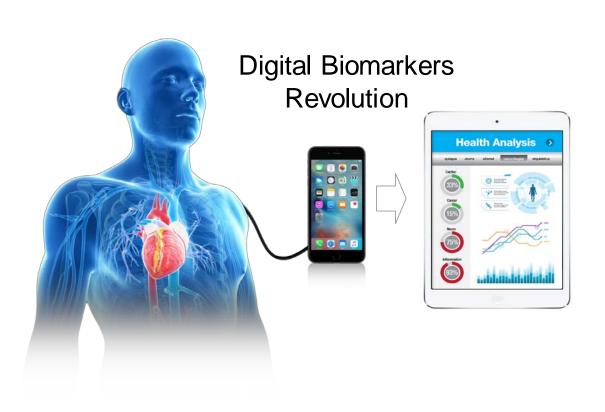
Q1 2020 Growth Stratification





Convergence of Technology and Healthcare





Late: Invasive











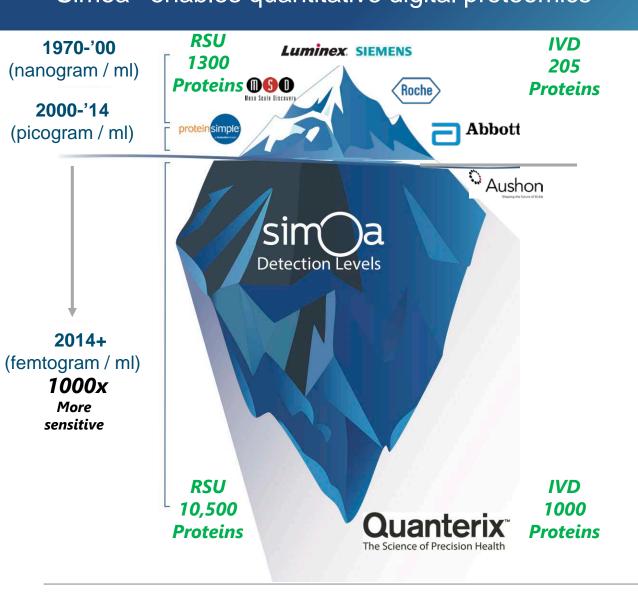


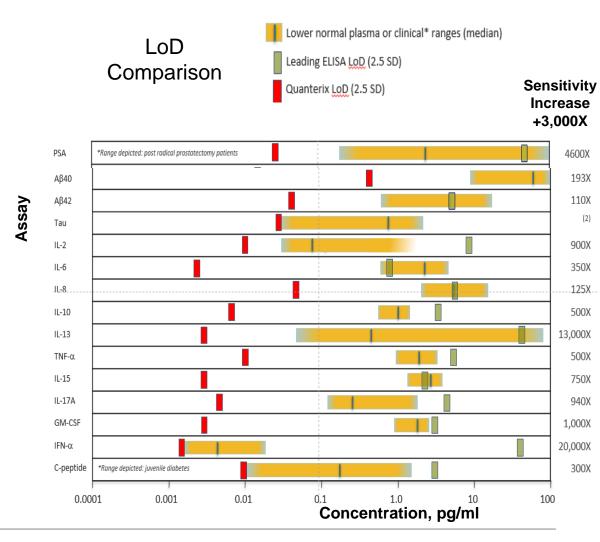


Early: Non-Invasive

Evolution of Immunoassay Simoa® enables quantitative digital proteomics







Compelling and Rare Investment Thesis Digital Proteomics Disruption



EXECUTION

Low Risk Research
Short to Mid-term: 2 – 3x

"High Probability" Drug Development

- ✓ Strong cash position; multiple growth catalysts
- ✓ Global PPH ecosystem engaged for COVID-19
- √ Validated; low regulatory/reimbursement risk
- ✓ Strong economic profile, early penetration
- ✓ Rapid Adoption \$0 to \$57M as of 2019
- ✓ Proven mgt, board and founder

ASPIRATION

High Value Diagnostics & Health Screens
Mid to Long-term: 10 – 15x

"Liquid MRi" for Brain, Serology & "Liquid Biopsy"

- ✓ Own100% rights; CDx → LDT → IVD
- ✓ Disruptive blood test with high optionality
- ✓ Covid-19, Neuro, liquid biopsy, drug rescue
- ✓ High Fidelity Serology test under consideration
- ✓ High potential for partnerships in COVID / Neuro
- ✓ Abbott, Quest, Labcorp, Siemens, Roche, Myriad

Capture Major Market Opportunity





\$1B - \$4B: Research

Market Growth ~10% p.a.

Oncology / Immunology

Luminex (55% Share) MSD (17%) Others (28%)

HD-X & SP-X < 1% Penetrated

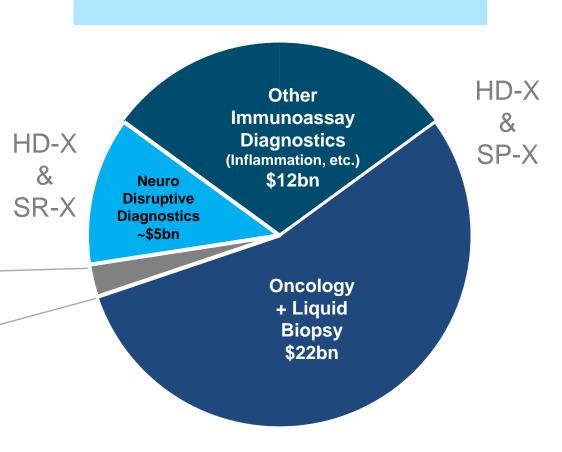
Neurology Low Competition

HD-X & SR-X
< 10%
Penetrated

Neuro
\$350m
Onco
\$750m

Body of evidence

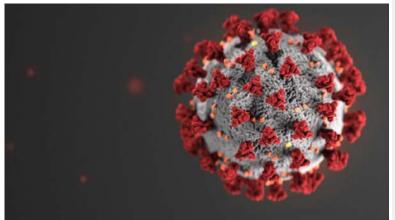
ASPIRATION \$40B+ Diagnostics



Neuro + Near-term CAC PIVOT



COVID-19 / ID



Catalyzed PPH global ecosystem:

- Measure total immune response
- Drug Targets and study Cytokine Storm
- **Simoa High Definition Serology RUO**

Accelerator Services



Expanded capacity and capabilities of Accelerator Laboratory services to:

- Overcome customer interruptions
- Sustain clinical trials given breakthroughs utilizing biomarkers drug approvals
- COVID emergency testing needs

China



Accelerate growth as people return to work:

- Drive rapid adoption in COVID-19 Neurology, Oncology, Inflammation
- Build local Accelerator Services Lab

COVID-19 Active Collaborations PPH Network enabling worldwide outreach





COVID-19 High Resolution Test Suite: The Complete Picture Catalyzing PPH Ecosystem to Accelerate Commercialization





Innate Immune Response

- Cyto/Chemokines; HDX, SRX, & SPX
- Cytokine storm, vaccine & drug dev'l



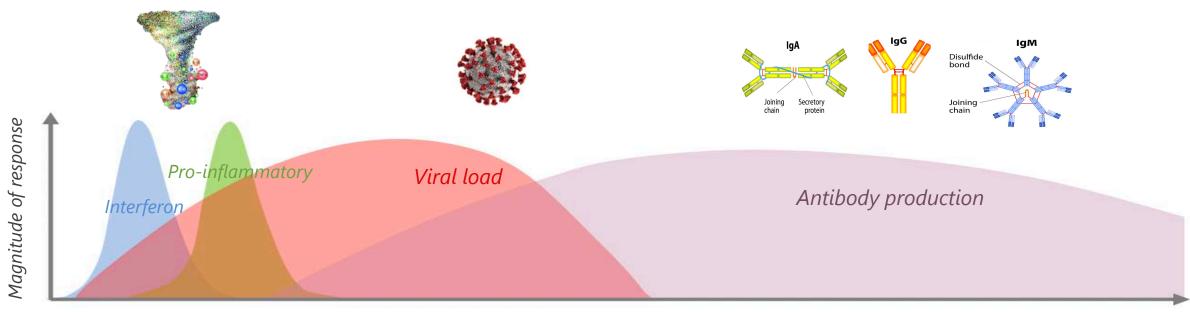
Viral Load

- Blood & saliva; HDX & SPX
- Apps: RUO & EUA for drug dev'l



Adaptive Immune Response – Serology

- High Def; HDX: EUA & HDX / SPX: RUO
- Immunity Passport, Disease severity, vaccine & drug dev'l, convalescent serum



Measure Innate Immunity Dysfunction Cytokine Storm radar system



Ultra sensitive cytokine & chemokine panels to help anticipate & manage immune dysregulation

CorPLexTM Cytokine Panel: IFN-γ, IL-1β, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-22, TNFα

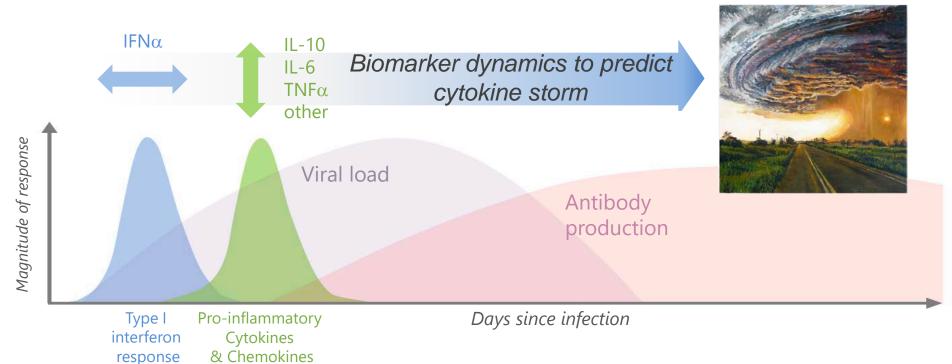
Chemokine Panel IP-10, ITAC, MCP-1, MIP-3b

Cytokine 6-Plex Panel IFN-γ, IL-6, IL-10, IL-12p70, IL-17A, TNFα



Simoa HD-X

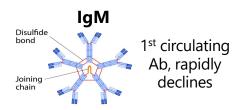
Simoa SP-X



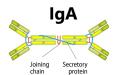
- Predict disease severity and stratify patients into risk groups
- Administer tailored immune modulation therapy and monitor response
- Vaccine, Drug and Convalescence plasma treatment development

QTRX COVID-19 High Definition Serology Panel





Most abundant, long-term immunity



Mucosal immunity, pathogen clearance



- Comprehensive immunity: IgG 1st & can add IgM / IgA
- First quantitative IgG test in serum/plasma in development
- High sensitivity / specificity; Reduce false +/-'s, See early
- Home (finger prick, dry blood spots) being explored
- High throughput (1000 tests/inst/day) for large scale testing
- Global population testing via PPH, instr's and sample banks
- Validation underway; Goal RUO launch date: IgG Q2'20

Immune response

Susceptibility and recurrence risk

Early viral detection

Viral carriers/ asymptomatic

Active Immunity

Passive immunity and vaccine development

Viral evolution, future protection



Innate Immune Response

• Simoa COVID-19 Cytokine & Chemokine panels for cytokine storm signatures, vaccine & drug development



Viral Load

Direct viral detection in **blood** & saliva to support diagnosis & drug development



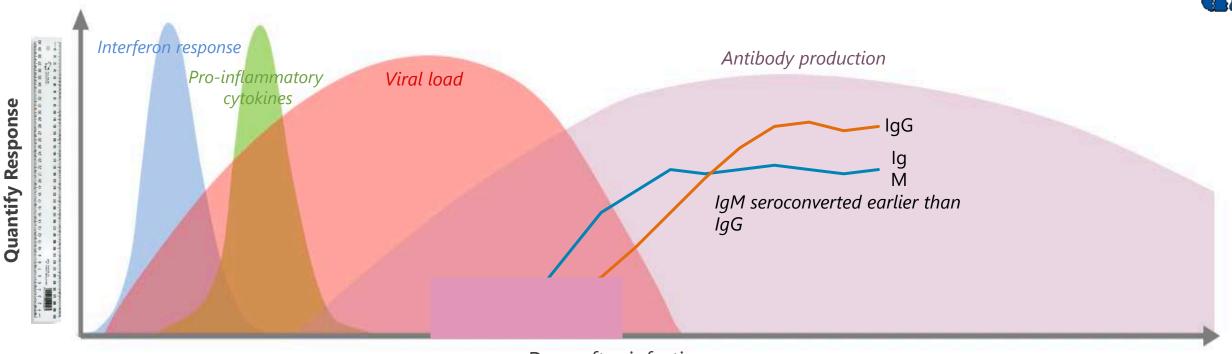
Adaptive Immune Response Profiling

- Simoa COVID-19 lgG & High Definition **Serology** for Immunity, Severity, vaccine & drug dev'l, convalescent serum qualification
- Low-invasive, quantification of seroconversion













Innate Immune Response

• Simoa COVID-19 Cytokine & Chemokine panels for cytokine storm signatures, vaccine & drug development



Viral Load

Direct viral detection in blood & saliva to support diagnosis & drug development



Adaptive Immune Response Profiling

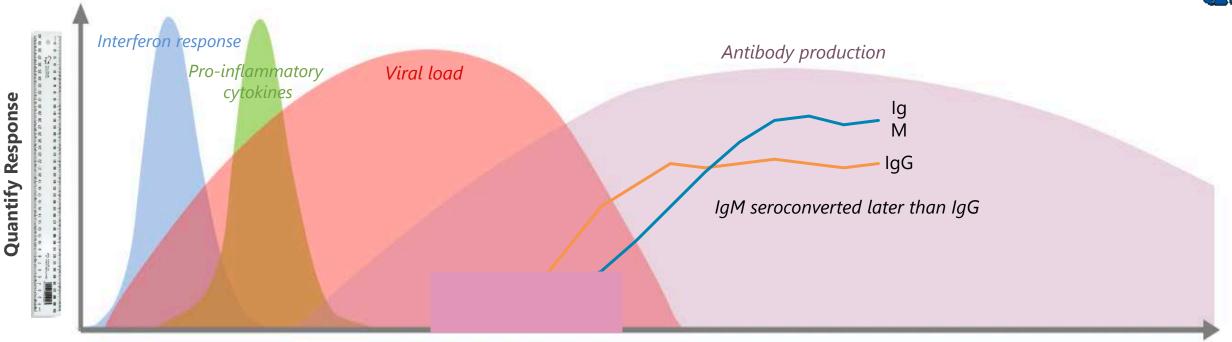
- Simoa COVID-19 lgG & High Definition **Serology** for Immunity, Severity, vaccine & drug dev'l, convalescent serum qualification
- Low-invasive, quantification of seroconversion







1. Quantitation







Quantify Response

• Simoa COVID-19 Cytokine & Chemokine panels for cytokine storm signatures, vaccine & drug development





Direct viral detection in **blood** & saliva to support diagnosis & drug development

Adaptive Immune Response Profiling



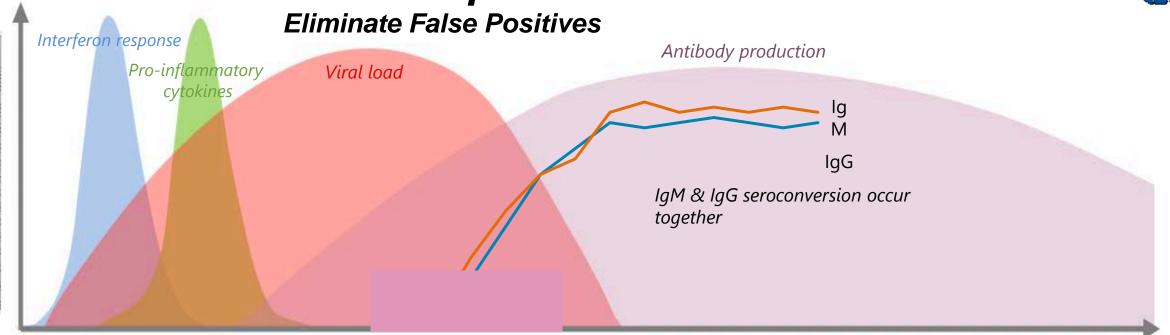
• Low-invasive, quantification of seroconversion







2. Ultra Specific







Quantify Response

Innate Immune Response

Simoa COVID-19 Cytokine & Chemokine panels for cytokine storm signatures, vaccine & drug development



Viral Load

Direct viral detection in blood & saliva to support diagnosis & drug development



Adaptive Immune Response Profiling

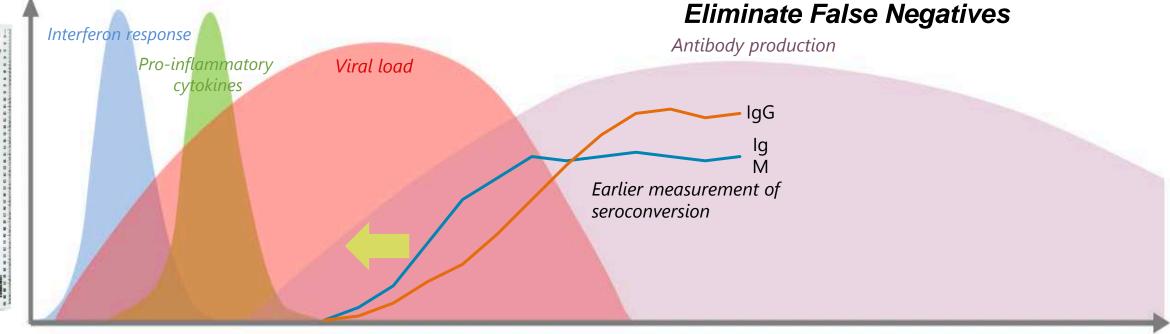
• Simoa COVID-19 lgG & High Definition **Serology** for Immunity, Severity, vaccine & drug dev'l, convalescent serum qualification







3. UtraerSensitive





1

Quantify Response

Innate Immune Response

 Simoa Cytokine & Chemokine panels for cytokine storm signatures, vaccine & drug development



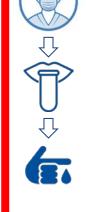
Viral Load

Direct viral detection in blood & saliva to support diagnosis & drug development

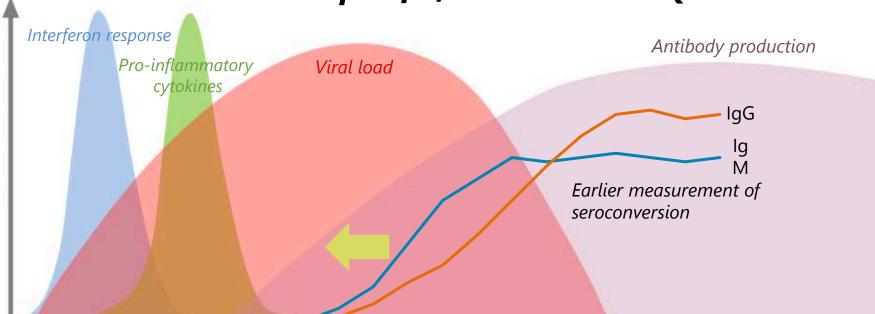


Adaptive Immune Response Profiling

• Simoa lgG & **High Definition Ruo Serology** for Immunity, Severity, vaccine & drug dev'l, convalescent serum qualification



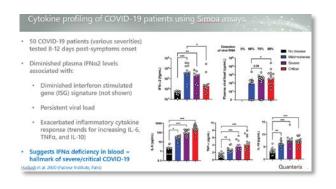
Ultra Specific, Sensitive and Quantitative



Commercial CAC - Pivot Plan COVID19 Immune Response, Drive Accelerator and Expand China footprint



Program	Pivot Plan			
<u>C</u> OVID	 PPH Direct ID customer outreach in hot zones (42 meetings in April) Optimized lead gen process and leverage PPH webinar catalyst Cytokine storm Radar and drugs, vaccine and plasma dev'l 			
<u>A</u> ccelerator	 Active engagement in global COVID clinical trials – Vaccines Recruit immune response population, patient & drug studies Bus Dev / Product linkage + Neuro retrospective trial focus 			
<u>C</u> hina	 Expand sales, ID KOL outreach, webinars, install base utilization COVID Collaborations w/ top China universities and assay innovators Expand APAC distributor and PPH reach via China hub and app center 			





















Quanterix Product Offerings



Instruments

Assay kits

Services

Bead based









HD-X

List Price: \$200k



255+ Installed @ YE 2019 (200+ HD-1s, 55+ HD-Xs)

Fully automated

1-6 Plex

Neuro Focused



SR-X

120+ Installed @ YE 2019

Benchtop

1-6 Plex

Neuro Focused



SP-X

List Price: \$75k

15+ Installed @ YE 2019

Benchtop

1 - 10 + plex

Onco Focused



Plate

Bead

300+ assays

Homebrew kits

Singleplex and Multiplex

100 Plate Kits 80 Bead Kits



Accelerator

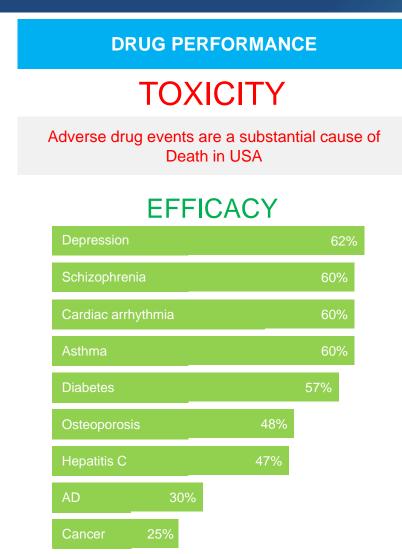
Contract research

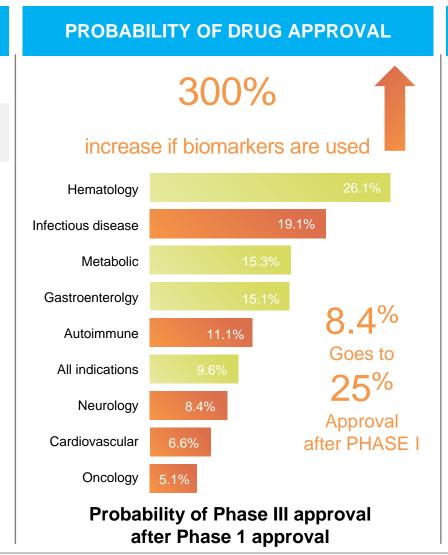
Custom assay development & reagent production

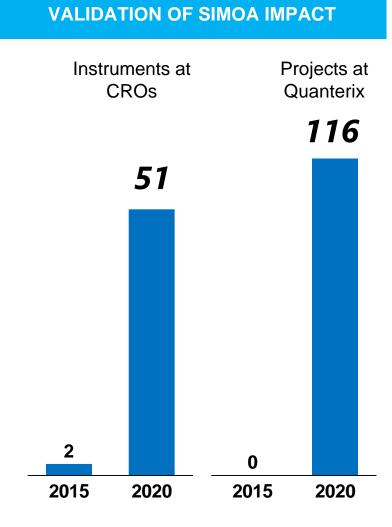
CLIA and **LDT** capabilities

Addressing a Significant Unmet Need in Drug Development



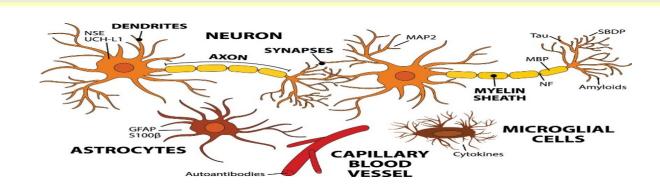






FDA Supporting Biomarkers for Drug Development





Neuro health is national Priority:

- veteran PTSD
- opioid addiction
- rapid AD ramp
- mental issues
- healthcare cost burden

Drug makers discouraged by lack of returns and subjective cognitive endpoints Issued guidance to use biomarkers for drug approvals

Rapid adoption

Source: Health Advances analysis, Quanterix materials.

Blue Chip Customers 2019



Research Institutions



VAN ANDEL

MGH

WAYNE STATE

Biopharma



CROs & Other























immun













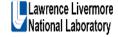




































RESEARCH







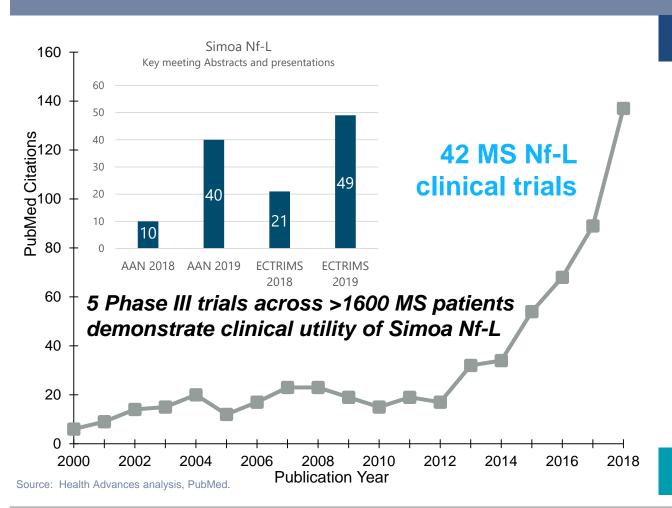




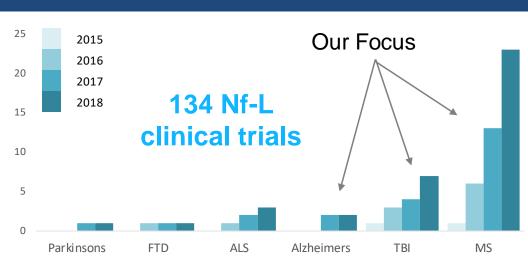
Nf-L Rapidly Expanding as Best In Class Neuro Biomarker



PEER-REVIEWED PUBLICATIONS ON Nf-L



Nf-L PUBLICATIONS



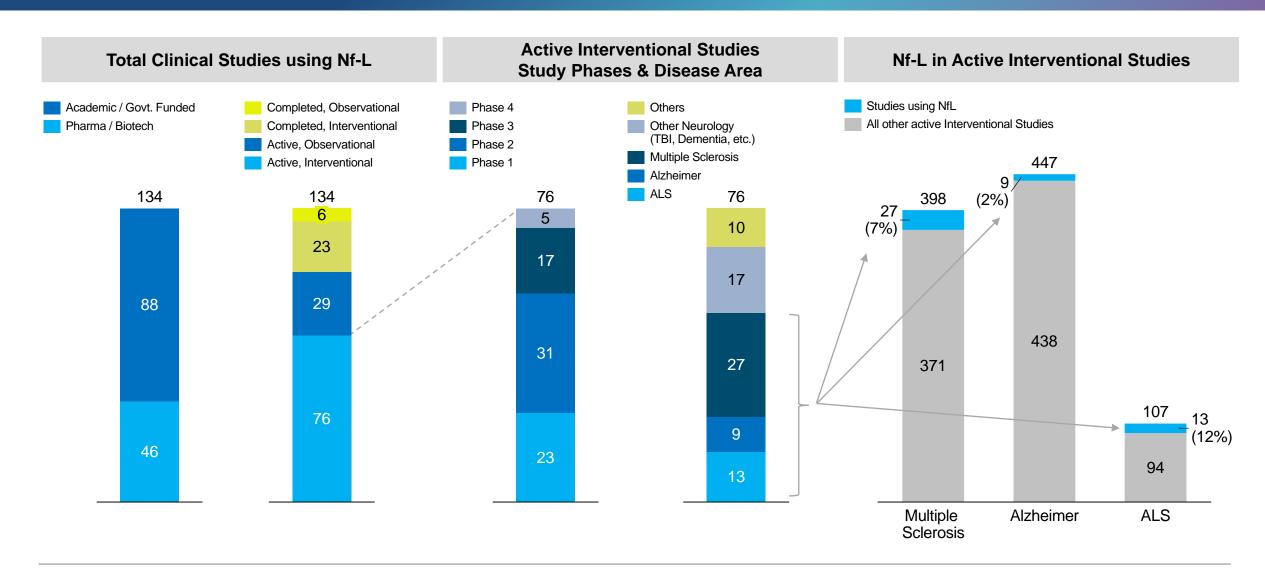
STUDIES CONFIRM NFL CLINICAL UTILITY:

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

Majority of published data obtained with Simoa Nf-L

Unlocking rich clinical potential of Nf-L





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













No disability with only

minimal signs

Moderate disability

Disability affects full daily activities Essentially restricted to wheelchair

Drug Trials





- Phase III ASCLEPIOS I/II trials
- Nf-L used as secondary endpoint, reduction after 3 months

Drug Selection





- Phase III OPERA I/II AND ORATORIO trials: **Nf-L lowered to healthy.**
- Ocrevus: \$1.75B sales, highest MS drug share in US of 40%

Disease Biology



- Understanding of MS genetics
- sNf-L increased 6 years before onset of MS

Advancing Toward Clinic

JAMA Neurology



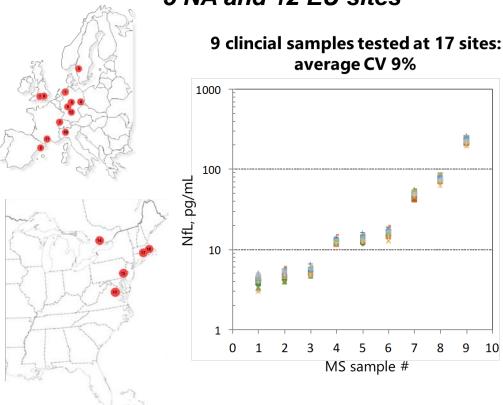


 sNf-L potential IVD assay, drug selection, drug / disease monitoring and health screens.

Neurofilament light chain (Nf-L): From Promising Biomarker to Clinically-relevant Tool

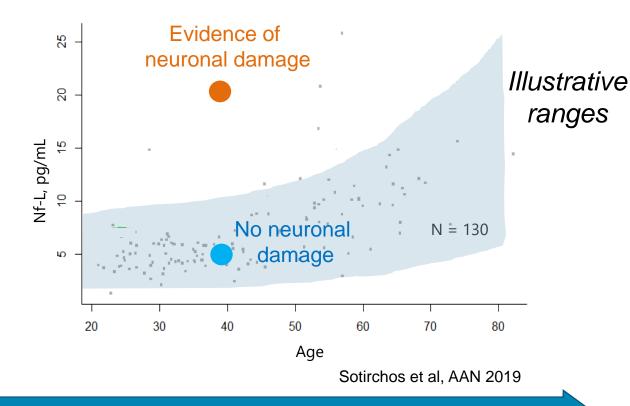






Large normative study ongoing 11,000 healthy controls, >18,000 samples





Analytical validity

Clinical Validity

2020 Objectives Revised in-light of COVID-19

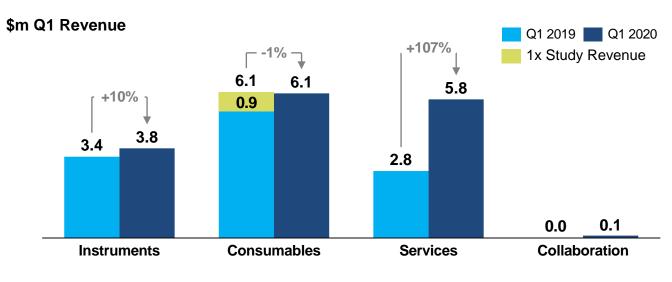


	Neurology	COVID-19 / ID	Strategy	Financials	Technology
40%	High double digit growth 40% 50%+ utilization	Develop Solutions for High-Resolution COVID- 19 Immune Response \$3-5m Revenue from JOLT-PIVOT	IVD partnerships Strategic M&A Diagnostics Strategy	3yr Growth: 30-40% Gross Margin: +200 bps Instr Rev Growth: 25%+ 75 90 HD-X's 75 90 SR-X/SP-X	100x sensitivity increase by YE 2021

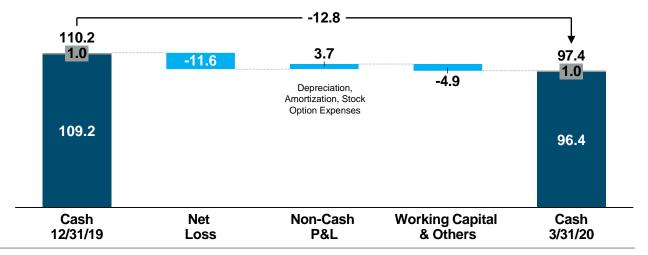
Q1 2020 Financial



In \$m	GAAP		Non-GAAP*		
	Q1 2020	Q1 2019	Q1 2020	Q1 2019	
Instruments	3.8	3.4	3.8	3.4	
Growth vs. PYR	10%		10%		
Consumables	6.1	6.1	6.1	6.1	
Growth vs. PYR	-1%		-1%		
Services	5.8	2.8	5.8	2.8	
Growth vs. PYR	107%		107%		
Collaboration	0.1	0.0	0.1	0.0	
Total Revenue	15.7	12.3	15.7	12.3	
Growth vs. PYR	27%		27%		
Gross Profit	6.8	6.0	7.6	6.0	
Gross Margin %	43.3%	48.7%	48.5%	48.7%	
Operating Expenses	18.5	15.4	18.4	15.2	
Loss from Operations	(11.7)	(9.4)	(10.8)	(9.2)	



\$m YTD Cash Flow



^{*} Non-GAAP item. Reconciliations are included in the Appendix to this presentation.

Poised to Disrupt Healthcare and Create Significant Value



Value

Execution: 2 – 3x value creation

Aspirational: 10 – 15x value creation

Best in Category-defining; Unrivaled Sensitivity / Technology Disrupt Class \$3B to New Methodical market penetration strategy to reward investors \$40B Answers market DNA – RNA - Protein; Better linked to Disease / Health Holy Grail Quanterix 700+ pubs Proven Validation: 19/20 top pharma, PPH, Drug Trials All Areas

penetration

5 Growth & Value; Razor – razor blade, \$150M invested

6 Low Risk / Solid Return + Uber Return Prospect

7 Track Record for Commercializing Disruption



Product

Launches

Differentiator

Rapid

Growth

Retail

Lynchpin

Quanterx

APPENDIX

Reconciliation of non-GAAP Financials in thousands (unaudited)



	2020		2019	
	Three months ended			
	Mar	ch 31		
Total revenue	\$ 15,727	\$	12,337	
GAAP Gross profit	\$ 6,813	\$	6,007	
Acquisition-related purchase accounting charges (Note 1)	\$ 820		_	
Non-GAAP gross profit	\$ 7,633	\$	6,007	
GAAP gross margin %	43.3%		48.7%	
Non-GAAP gross margin %	48.5%		48.7%	
GAAP total operating expenses	\$ 18,541	\$	15,364	
Acquisition-related purchase accounting charges (Note 2)	\$ 150	\$	151	
Non-GAAP total operating expenses	\$ 18,391	\$	15,214	
GAAP loss from operations	\$ (11,728)	\$	(9,357)	
Non-GAAP loss from operations	\$ (10,758)	\$	(9,207)	

Note 1: During the three months ended March 31, 2020, we incurred \$438 thousand of acquisition-related amortization of inventory valuation in connection with our acquisition of UmanDiagnostics. Also during the three months ended March 31, 2020, we incurred \$382 thousand of acquisition-related amortization of intangible assets in connection with our acquisition of UmanDiagnostics.

Note 2: During the three months ended March 31, 2020, we incurred \$150 thousand of acquisition-related amortization of intangible assets in connection with our acquisitions of UmanDiagnostics and Aushon Biosystems. During the three months ended March 31, 2019, we incurred \$151 thousand of acquisition-related amortization of intangible assets in connection with our acquisition of Aushon Biosystems.