

Q3 2021 Earnings Call

Kevin Hrusovsky, Chairman and CEO November 4, 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. 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. 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 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



Q3 2021 Advances

- Accelerated Growth
- SIMOA enabling therapeutic drugs



Financial Results

- Q3

- Cash flow



Objectives 2021



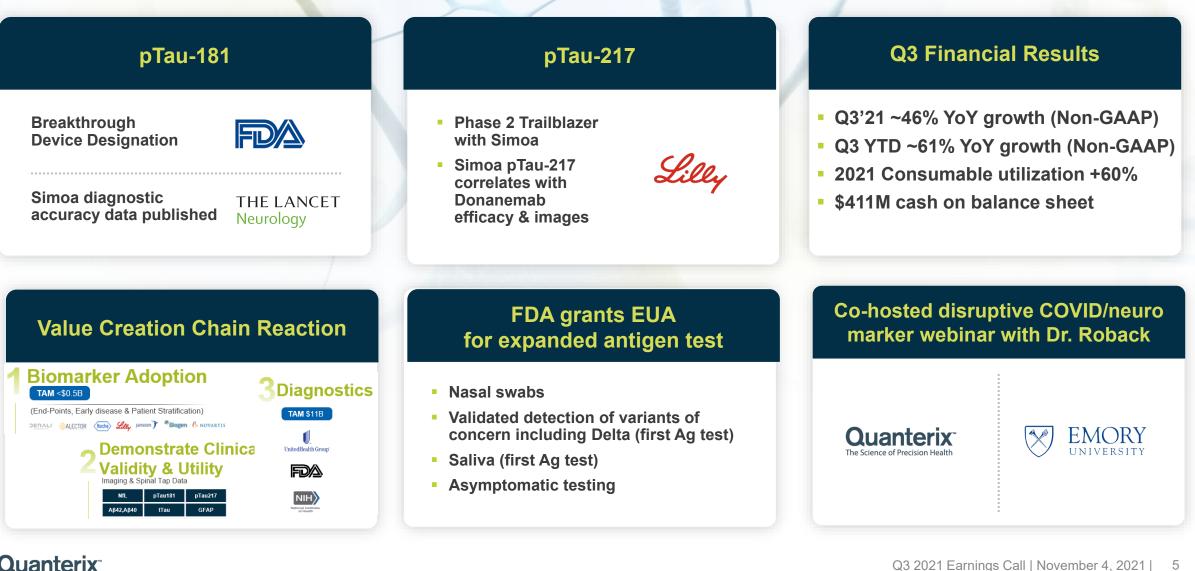
Blood-based

biomarkers

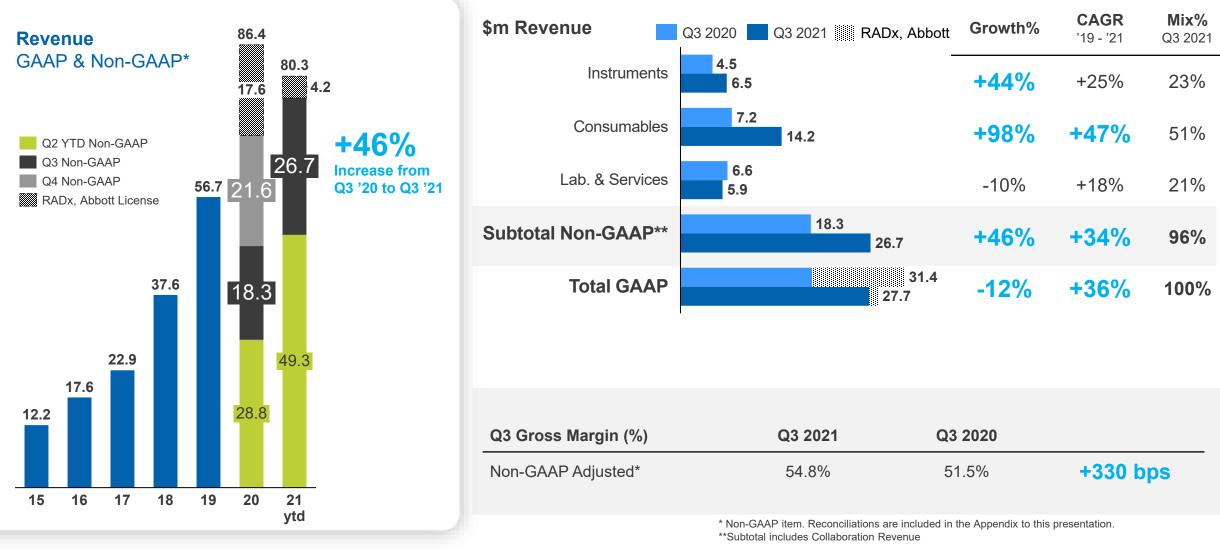
are fast becoming gold standard surrogates for traditional Tau or Abeta PET imaging Ever since the FDA approved Biogen and Eisai's Alzheimer's disease treatment Aduhelm, the buzzword among competing companies has been **"biomarker."** And that's just what Eli Lilly has to offer with two new analyses on their candidate donanemab."

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





Q3 2021 – Revenue Growth & Gross Margin

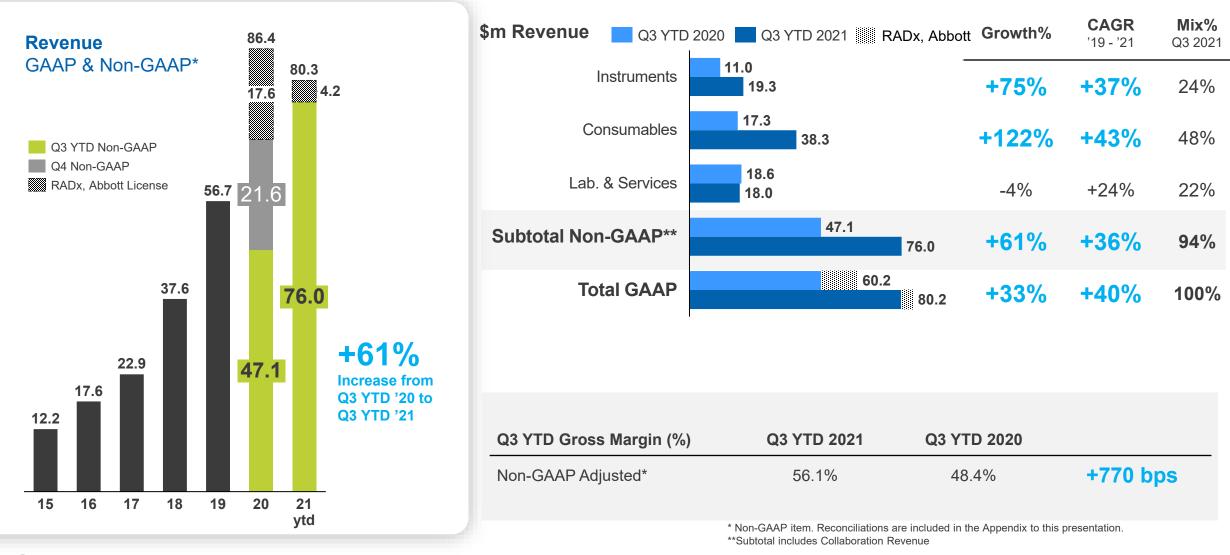


The Science of Precision Healt

6

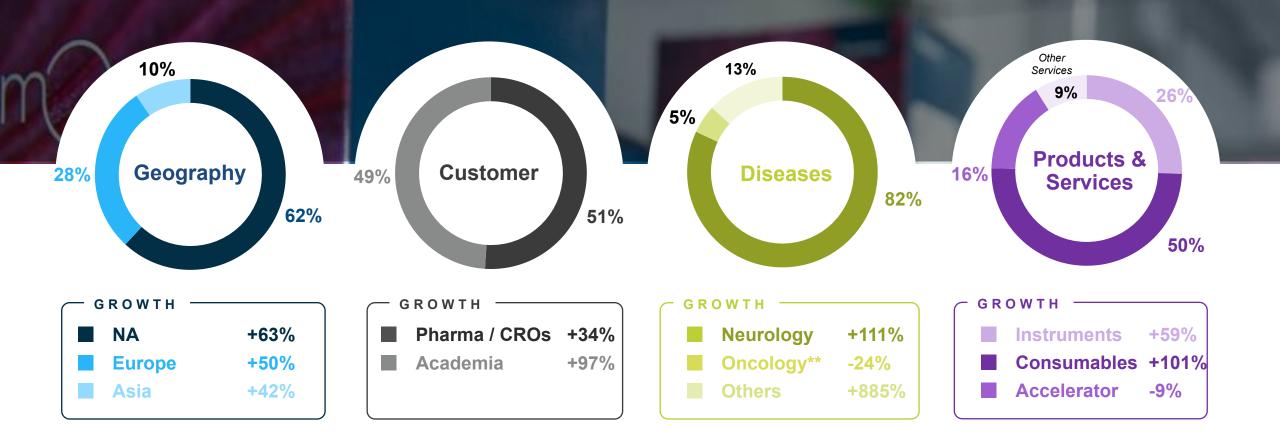
Q3 2021 ADVANCES

Q3 YTD 2021 - Revenue Growth & Gross Margin



Quanterix^{**}

Revenue Stratification (TTM)





Q3 2021 Advances Neurology Immunology & Oncology Infectious Diseases Others PUBLICATIONS **BIOMARKERS INSTRUMENTS** ACCELERATOR Cumulative Cumulative **Placements** Projects & Revenue 1,485 \$m Revenue # of units placed 18.0 Cumulative Number of Drug 1,120 Trial Projects 11.5 10.0 1.7 7.7 3.6 4.9 3.1 6.1 1.1 ytd ytd ytd ytd

Scientific Validation driving Adoption

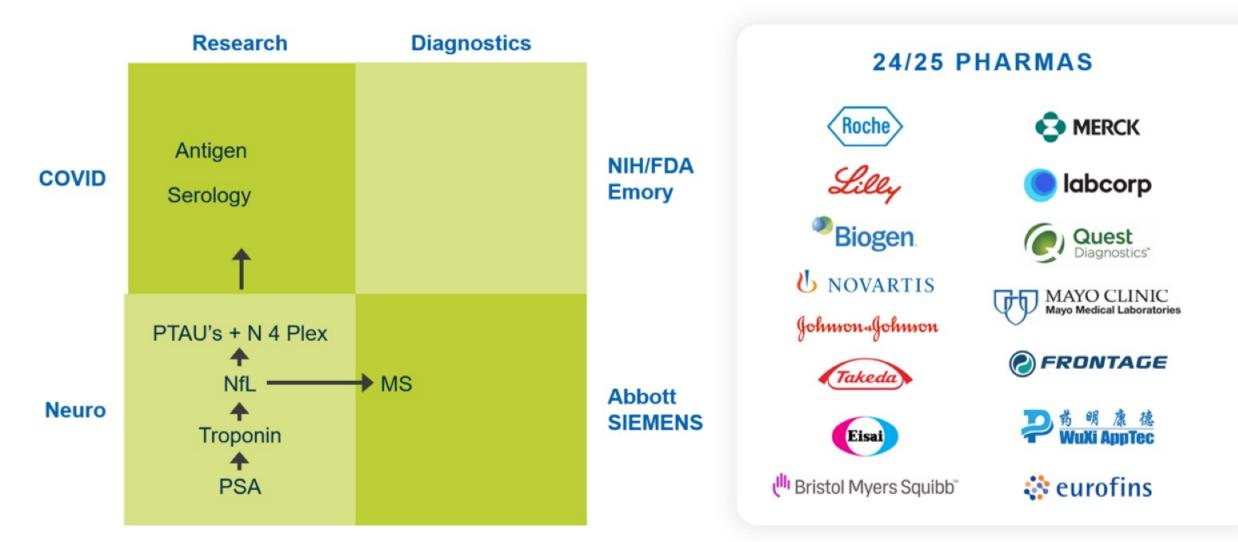
Q3 2021 Advances

Quanterix

The Science of Precision Heal

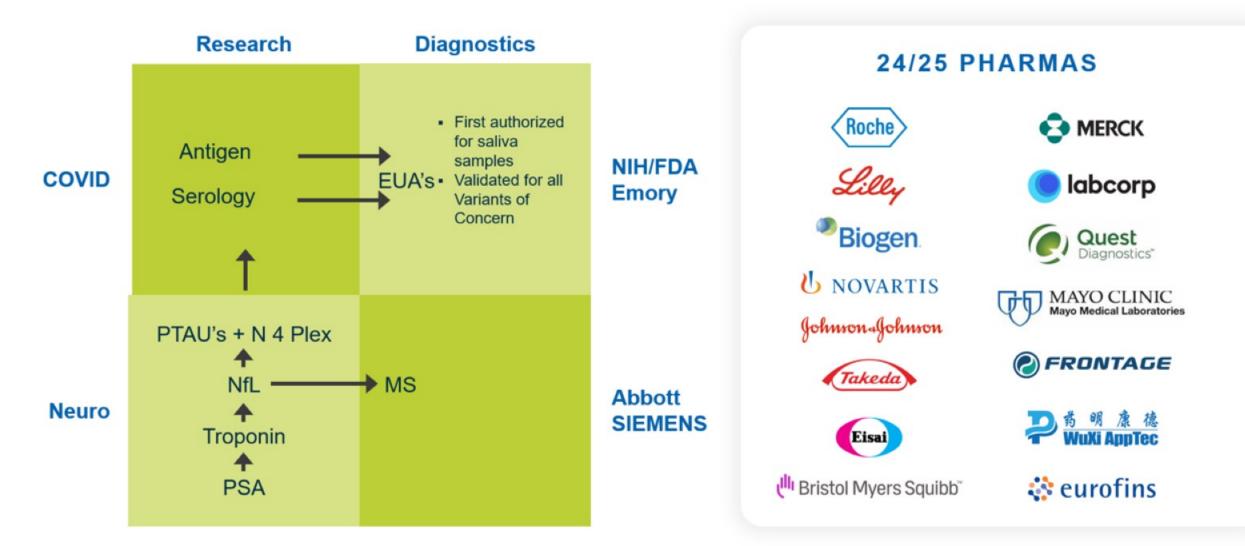
Q3 2021 Earnings Call | November 4, 2021 | 9

STRATEGIC ROAD MAP

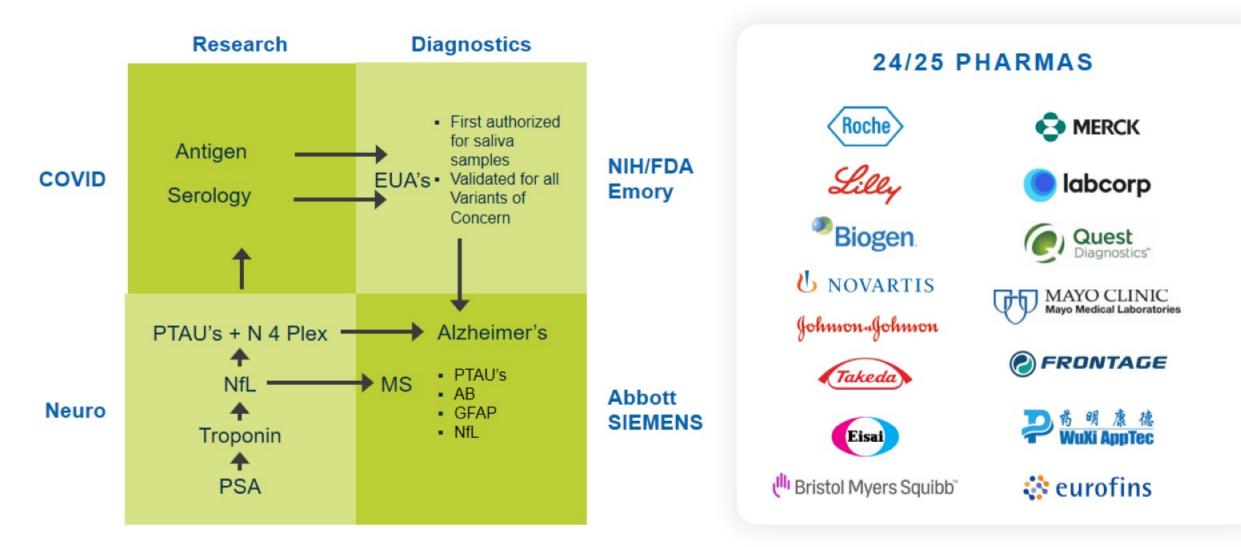




STRATEGIC ROAD MAP



STRATEGIC ROAD MAP





Compelling Thesis to Lead Proteomics into Precision Health Era

Execution

2 to 3x Value Creation

RESEARCH & DISCOVERY

\$1B to \$20B TAM²



- Strong Validation pTau181 Breakthrough Device Designation
- Strategy: Fortify Moat w/ 100x sensitivity for Neuro

Growth Catalysts

- pTau's & Neuro Plex: AD Diagnostic Therapies
- Revitalized Neuro drug trials
- Scale, HDx Clinical, Accelerator Expansion & 100x

Aspiration

10 to 15x Value Creation

DIAGNOSTICS & HEALTH SCREENS

Mounting Evidence – AAIC plasma pTau's

\$12B to \$100B TAM²

- 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

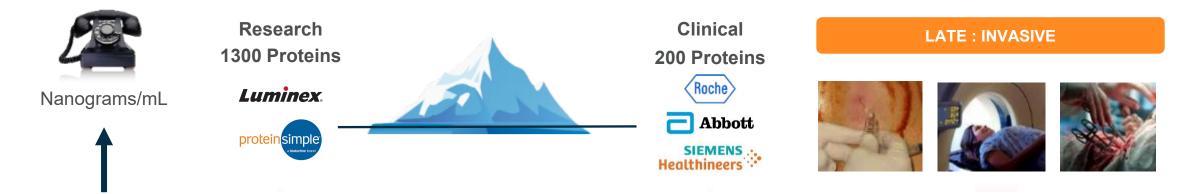
* 2017-2020 3Yr CAGR

1. 409A Valuation YE 2014

n YE 2014 2. Based on independent third-party research report 2021

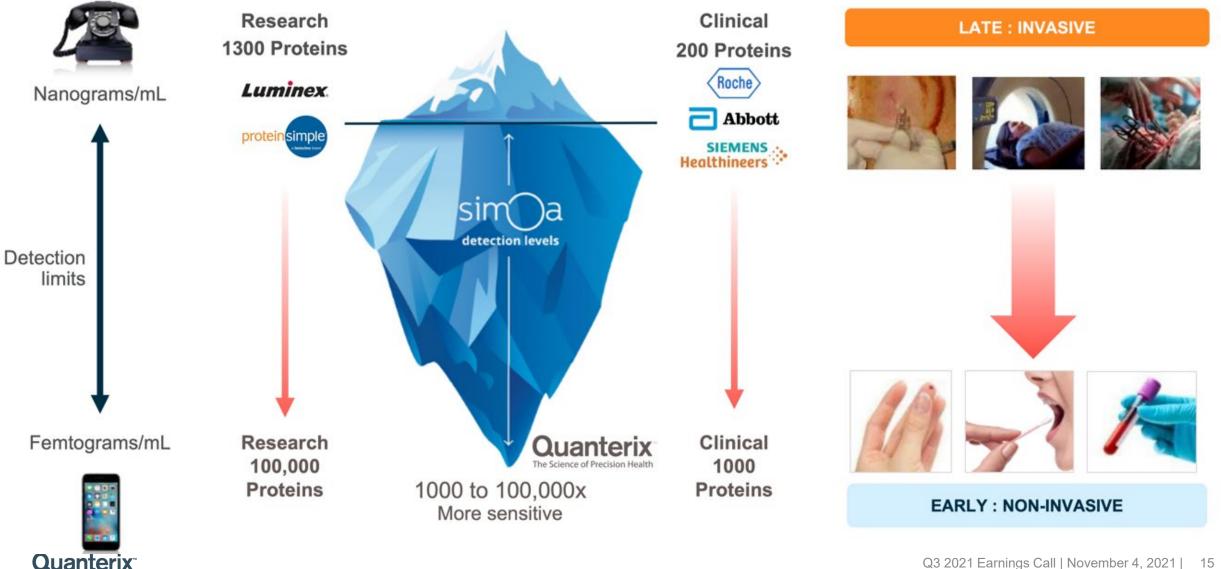


Digital Biomarkers Sensitivity Unlocking Proteomics

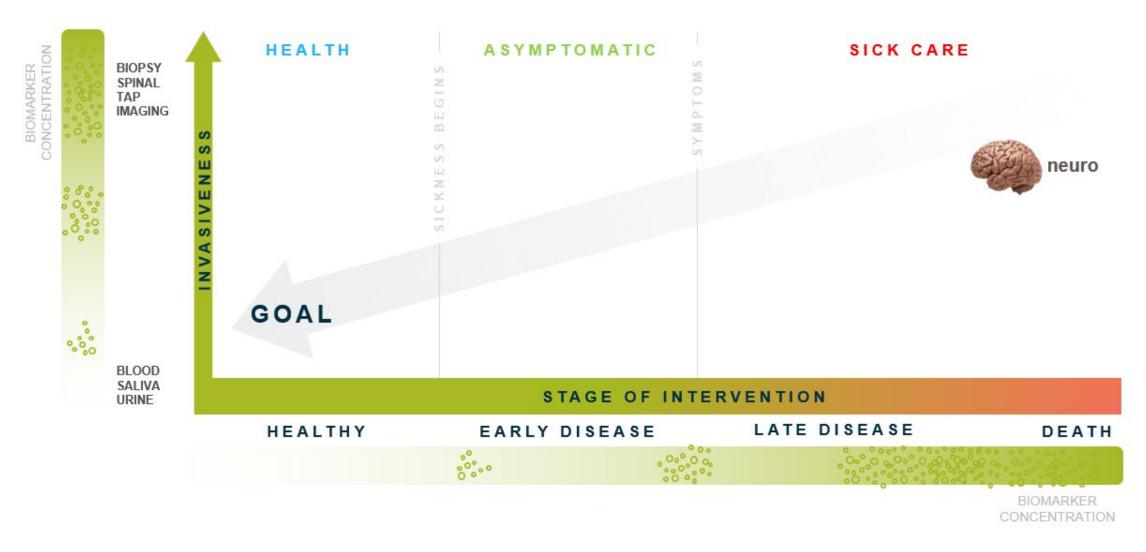




Digital Biomarkers Sensitivity Unlocking Proteomics

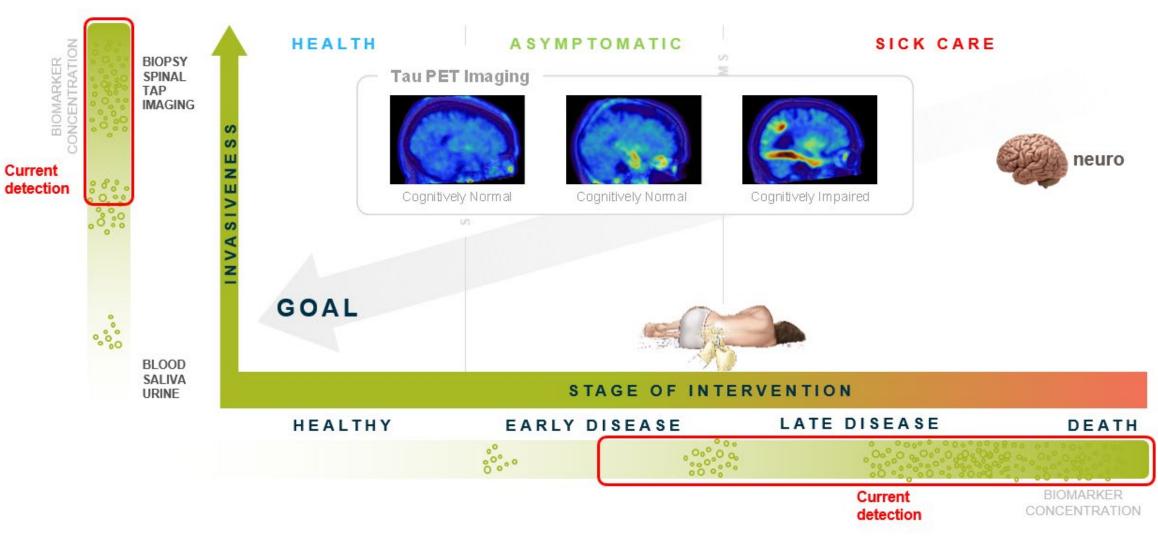


Simoa[®] provides insight into Health to Disease Continuum



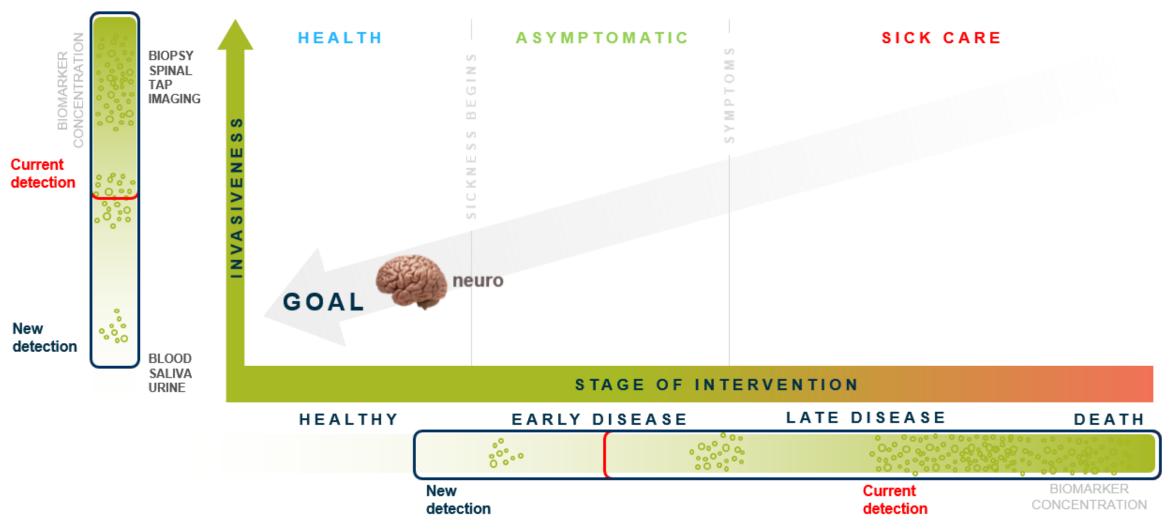


Simoa[®] provides insight into Health to Disease Continuum



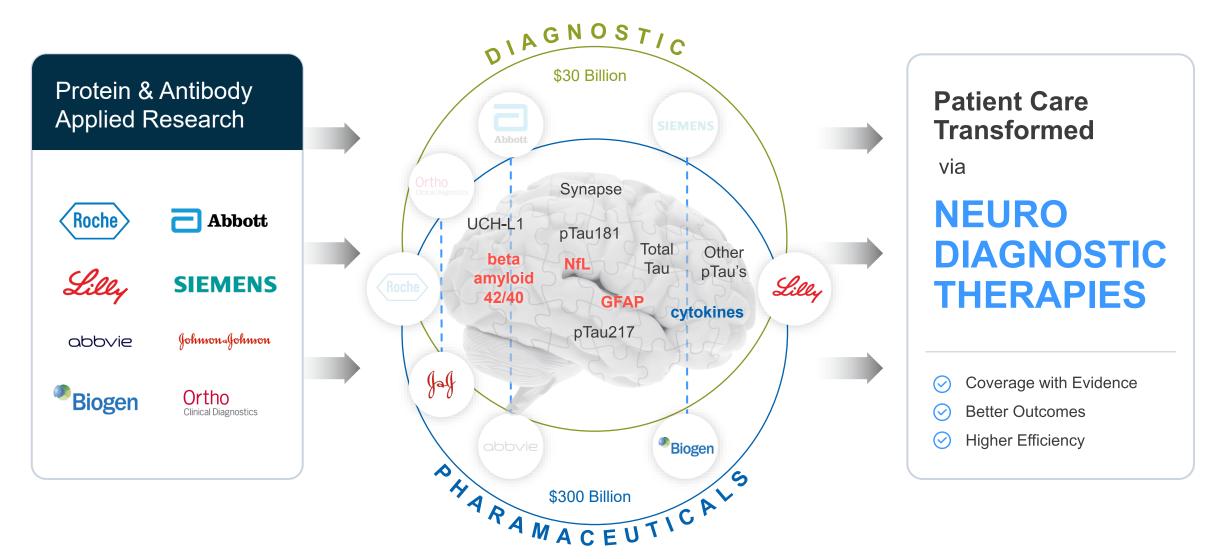


Simoa[®] provides insight into Health to Disease Continuum





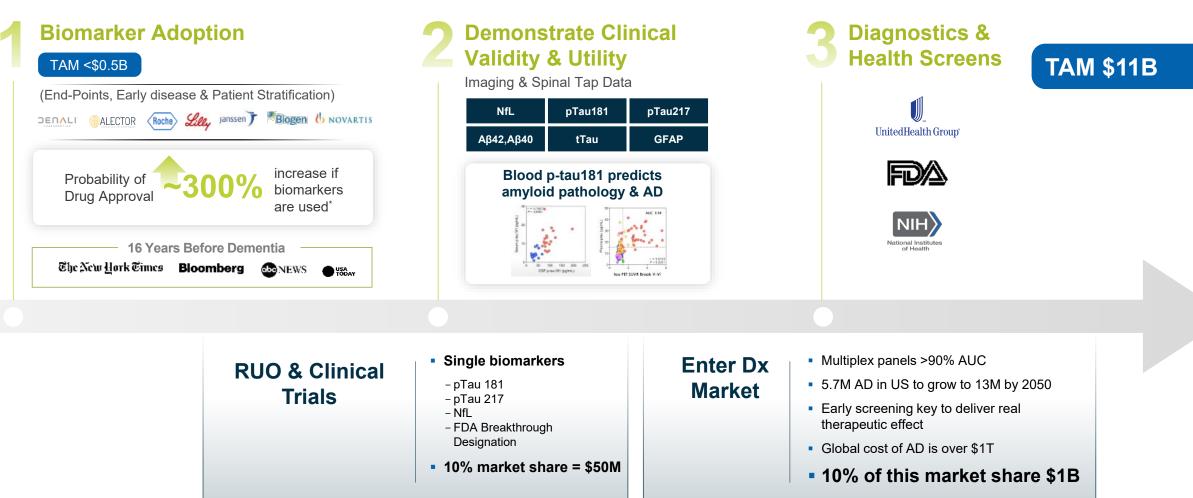
Potential for Establishing Category Through Diagnostic Disruption





Neurology poised for Value Creation Chain Reaction

Drug-trials center-piece of near-term focus



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





Eli Lilly Asks FDA to Approve Alzheimer's Drug

THE WALL STREET JOURNAL.

Home World U.S. Politics Economy Business Tech Markets Opinion Books & Arts Real Estate Life & Work WSJ. Magazine Sports ig Q

Eli Lilly Asks FDA to Approve Alzheimer's Drug

Drugmaker begins process seeking approval to sell donanemab as treatment for early-stage Alzheimer's



Tuesday, Lilly announces real-time submission for its donanemab to the U.S. FDA. Expects to complete its application in the next several months, which could lead to an FDA decision in the second half of 2022

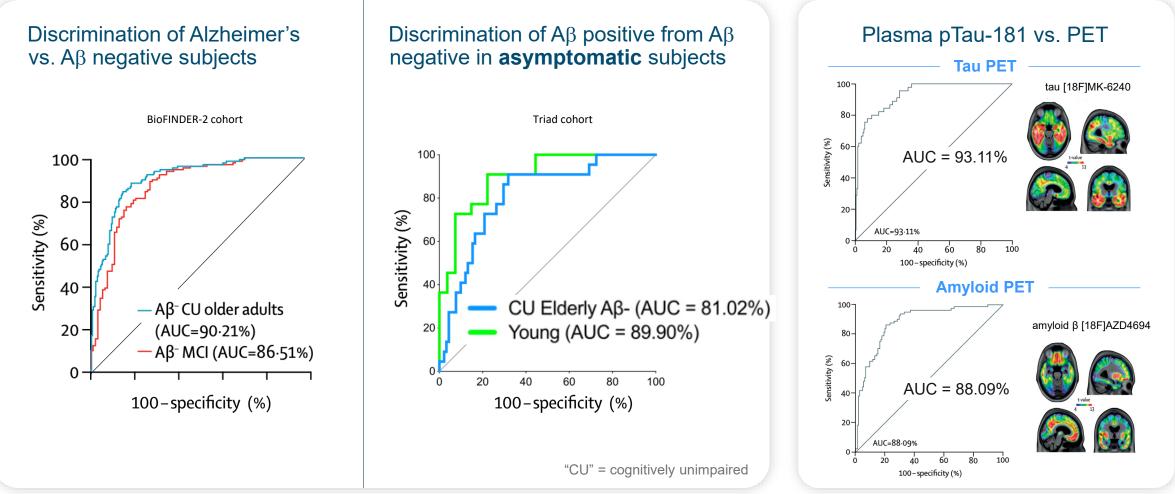
The Wall Street Journal, October 28, 2021



Source: https://www.wsj.com/articles/eli-lilly-asks-fda-to-approve-alzheimers-drug-11635262324

Simoa pTau-181 diagnostic accuracy

Representative breakthrough data published in Lancet Neurology¹

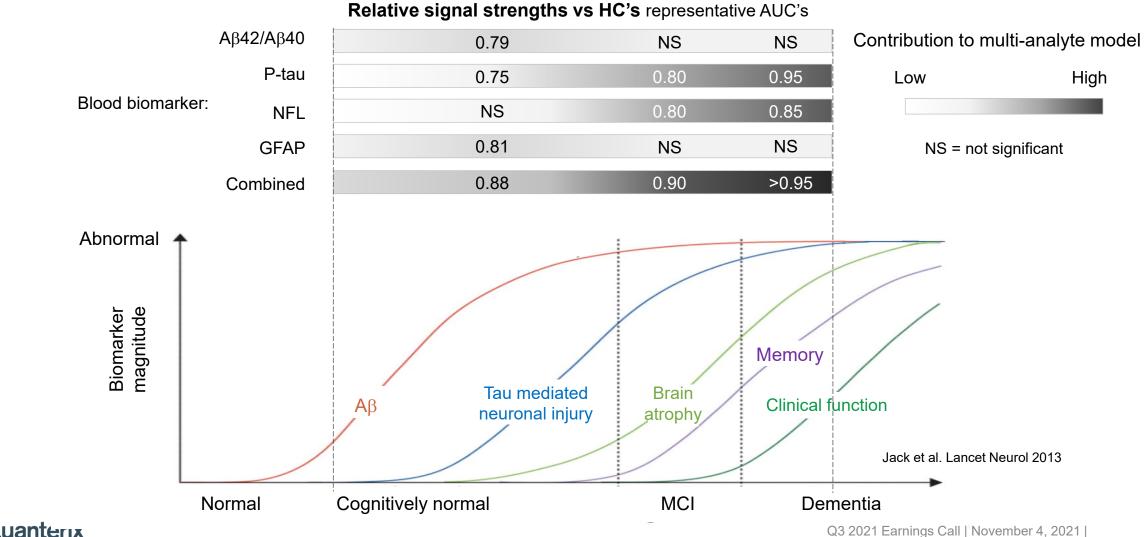


¹Karikari TK, Pascoal TA, Ashton NJ, et al. Lancet Neurol. 2020;19(5):422-433. doi:10.1016/S1474-4422(20)30071-5



Simoa multi-analyte approach to early Alzheimer's detection

Analyte signal contributions vary with the disease continuum



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



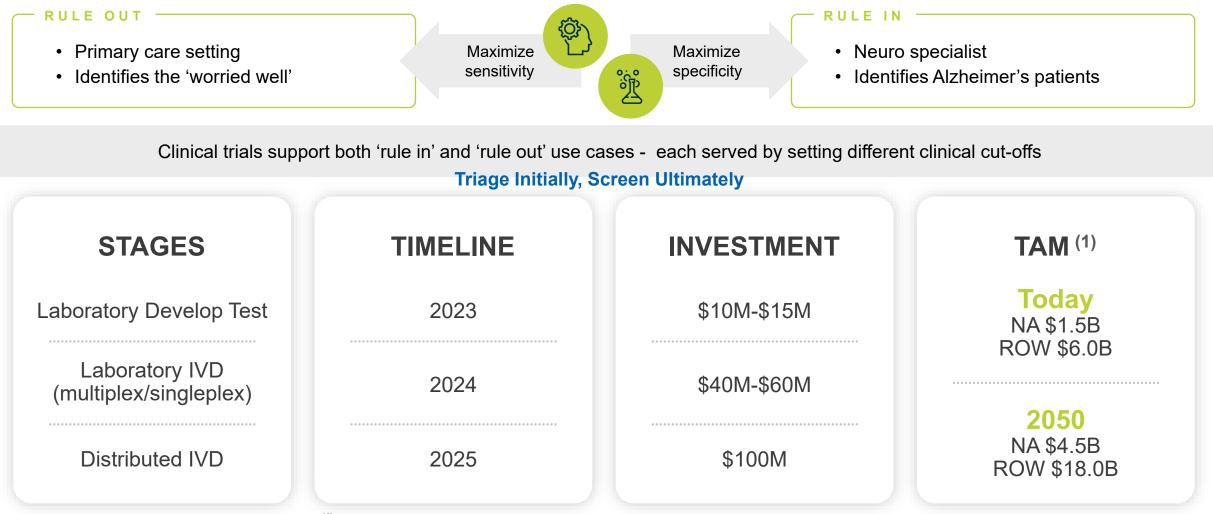
Source: www.alz.org



Accelerating Clinical Market Entry Alzheimer



Partnerships have the potential to accelerate and de-risk market penetration



⁽¹⁾ TAM assumptions are post dementia diagnosis of patients over 50 years of age with of 15 years of annual screening (20% NA / 80% ROW)

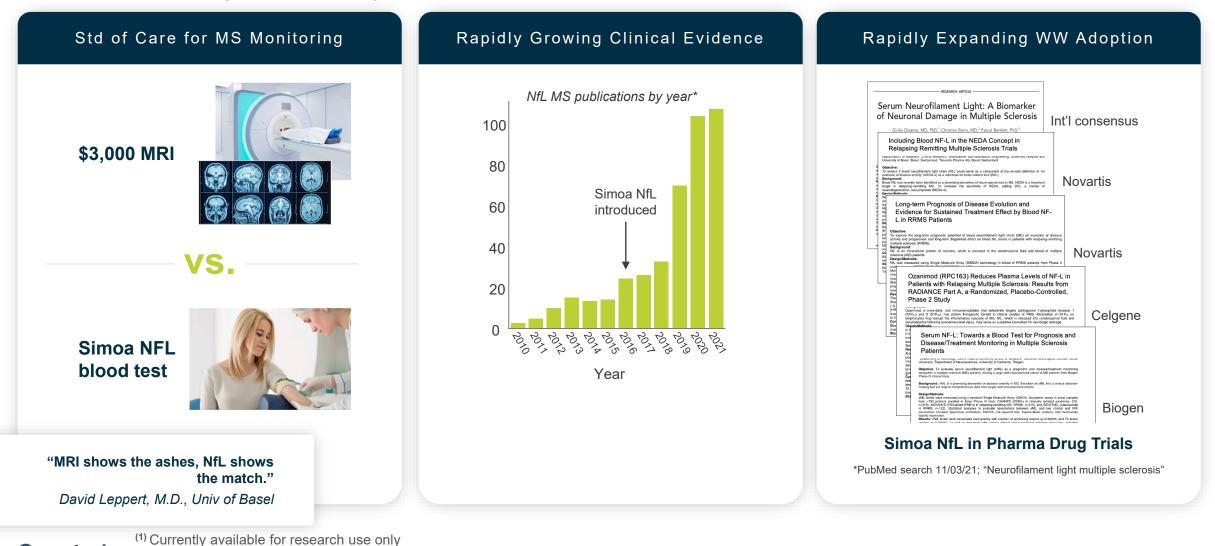
anterix

The Science of Precision Health

Source: PubMed

Simoa NfL in Multiple Sclerosis

QTRX has the only commercially available blood test for NfL⁽¹⁾



Q3 2021 Earnings Call | November 4, 2021 |

26

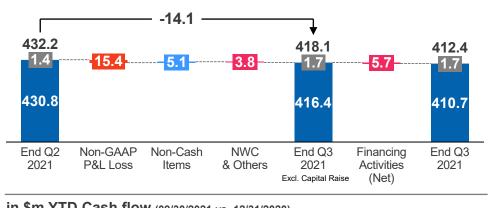
Q3 2021 Financials

	Q3 (3 Months)				YTD (9 Months)				
	GAAP		Non-GAAP*		GAAP		Non-GAAP*		
In \$m	2021	2020	2021	2020	2021	2020	2021	2020	
Instrument	6.5	4.5	6.5	4.5	19.3	11.0	19.3	11.0	
Growth vs. PYR	44%		44%		75%		75%		
Consumable	14.2	7.2	14.2	7.2	38.3	17.3	38.3	17.3	
Growth vs. PYR	98%		98%		122%		122%		
Product Revenue	20.7	11.7	20.7	11.7	57.6	28.3	57.6	28.3	
Growth vs. PYR	77%		77%		104%		104%		
Services	5.9	6.6	5.9	6.6	18.0	18.6	18.0	18.6	
Growth vs. PYR	-10%		-10%		-4%		-4%		
Collaboration & RADx	1.1	13.2	0.1	0.0	4.7	13.3	0.5	0.2	
Total Revenue	27.7	31.4	26.7	18.3	80.3	60.2	76.0	47.1	
Growth vs. PYR	-12%		46%		33%		61%		
Cost of Goods & Services	12.4	10.3	12.1	8.9	34.8	27.1	33.4	24.3	
Gross Profit	15.2	21.1	14.6	9.4	45.5	33.1	42.6	22.8	
Gross Margin %	55%	67%	55%	51%	57%	55%	56%	48%	
Operating Expenses	30.5	18.8	30.0	17.5	84.2	54.8	80.7	53.4	
Income/(Loss) from Operations	(15.2)	2.3	(15.4)	(8.1)	(38.7)	(21.7)	(38.1)	(30.6)	

- Record Product Revenue \$20.7m / +77%
- Non-GAAP Gross Margin 54.8%* /+330 bps vs. PYR, driven by volume, price and productivity

\$m Q3 Cash Flow

Restricted Cash



	Casil IIU	W (09/30/2021	vs. 12/31/202	20)	
182.6	-38.5	+13.8	-9.4	148.5	+264

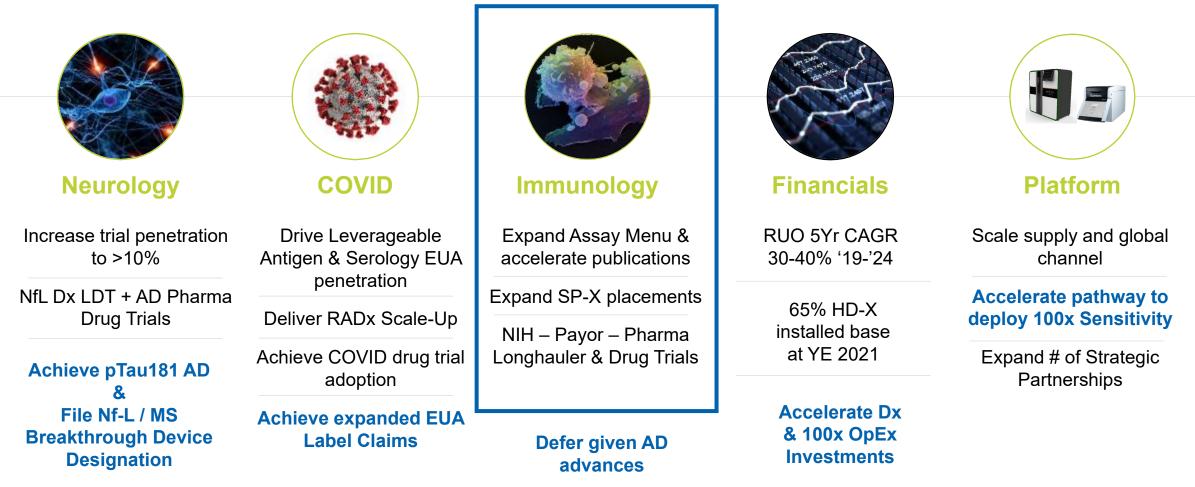
* Non-GAAP item. Reconciliations are included in the Appendix to this presentation.



412.4

Objectives 2021

RUO 2019-2024 CAGR 30-40%





APPENDIX





Use of Non-GAAP Numbers

	In \$m	Total Revenue	Cost of Goods Sold	Gross Profit	Gross Margin %	Operating Expenses	Loss from Operations
	GAAP	27.7	12.4	15.2	55.1%	30.5	-15.2
Q3 2021	Non-GAAP adjustments: Grant revenue (Note 1) Acquisition-related purchase accounting charges (Note 3)	-1.0	-0.4	-1.0 0.4		0.0	-1.0 0.4
	Grant research and development expenses (Note 5)					-0.5	0.5
	Non-GAAP	26.7	B Goods Sold Profit Margin % Experimentation 12.4 15.2 55.1% 30 -1.0 -1.0 -1.0 -0.4 0.4 -0.4 0.4 00 -0.6 0.4 12.1 14.6 54.8% 30 -0.6 34.8 45.5 56.6% 84 -1.4 1.4 -00 -3 33.4 42.6 56.1% 80 -1.9 -11.2 -0.4 0.4 00 -1.0 1.0 -0 -1 -1 8.9 9.4 51.5% 17 -1.8 1.8 -00 -1 -1.0 1.0 -0 -1	30.0	-15.4		
	GAAP	80.3	34.8	45.5	56.6%	84.2	-38.7
YTD 2021	Non-GAAP adjustments:						
	Grant revenue (Note 1)	-4.2		-4.2			-4.2
	Acquisition-related purchase accounting charges (Note 3)		-1.4	1.4		-0.1	1.5
	Grant research and development expenses (Note 5)					-3.4	3.4
	Non-GAAP	76.0	33.4	42.6	56.1%	80.7	-38.1
	GAAP	31.4	10.3	21.1	67.2%	18.8	2.3
	Non-GAAP adjustments:						
	Grant revenue (Note 1)	-1.9		-1.9			-1.9
	License agreement Revenue (Note 2)	-11.2		-11.2			-11.2
QJ 2020	Acquisition-related purchase accounting charges (Note 3)		-0.4	0.4		0.0	0.4
	Cost of license Revenue (Note 4)		-1.0	1.0			1.0
	Grant research and development expenses (Note 5)			0		-1.3	1.3
YTD 2021 Q3 2020	Non-GAAP	18.3	8.9	9.4	51.5%	17.5	-8.1
YTD 2020	GAAP	60.2	27.1	33.1	55.0%	54.8	-21.7
	Non-GAAP adjustments:						
	Grant revenue (Note 1)	-1.9					-1.9
	License agreement Revenue (Note 2)	-11.2		-11.2			-11.2
	Acquisition-related purchase accounting charges (Note 3)		-1.8			-0.1	1.9
	Cost of license Revenue (Note 4)		-1.0	1.0			1.0
	Grant research and development expenses (Note 5)					-1.3	1.3
	Non-GAAP	47.1	24.3	22.8	48.4%	53.4	-30.6

Note 1: During the three months ended September 30, 2021, we recognized \$1.0 million in revenue in connection with our workplan 2 award under the National Institute of Health Rapid Acceleration of Diagnostics Program. During the nine months ended September 30, 2021, we recognized \$4.2 million in revenue in connection with our workplan 2 award under the National Institute of Health Rapid Acceleration of Diagnostics Program. During the three and nine months ended September 30, 2020, we recognized \$1.9 million in revenue in connection with our workplan 1 award under the National Institute of Health Rapid Acceleration of Diagnostics Program.

Note 2: During the three and nine months ended September 30, 2020, we recognized \$10.0 million in license revenue in connection with a non-exclusive license agreement with Abbott Laboratories. Also, during the three and nine months ended September 30, 2020, we recognized \$1.2 million of previously deferred license revenue as a result of entering into the license agreement with Abbott Laboratories.

Note 3: During the three months ended September 30, 2021, we incurred \$382 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of UmanDiagnostics. During the nine months ended September 30, 2021, we incurred \$274 thousand of acquisition-related amortization of inventory valuation and \$1,148 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of umanDiagnostics, as well as \$60 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of UmanDiagnostics, as well as \$60 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of UmanDiagnostics recognized in operating expenses. During the three months ended September 30, 2020, we incurred \$40 thousand of acquisition-related amortization of inventory valuation and \$382 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of UmanDiagnostics. During the nine months ended September 30, 2020, we incurred \$671 thousand of acquisition-related amortization of inventory valuation and \$1,147 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of intangible assets adjustments in connection with our acquisition of intangible assets adjustments in connection with our acquisition of intangible assets adjustments in connection with our acquisition of intangible assets adjustments in connection with our acquisition-related amortization of inventory valuation and \$1,147 thousand of acquisition-related amortization of intangible assets adjustments in connection with our acquisition of umanDiagnostics recognized in operating expenses.

Note 4: During the three and nine months ended September 30, 2020, we incurred \$1.0 million in license fees in connection with our non-exclusive license agreement with Abbott Laboratories.

Note 5: During the three months ended September 30, 2021, we incurred \$461 thousand in research and development expenses in connection with our workplan 2 award under the National Institute of Health Rapid Acceleration of Diagnostics Program. During the nine months ended September 30, 2021, we incurred \$3.4 million in research and development expenses in connection with our workplan 2 award under the National Institute of Health Rapid Acceleration of Diagnostics Program. During the nine months ended September 30, 2021, we incurred \$3.4 million in research and development expenses in connection with our workplan 2 award under the National Institute of Health Rapid Acceleration of Diagnostics Program. During the three and nine months ended September 30, 2020, we incurred \$1.3 million in research and development expenses in connection with our workplan 1 award under the National Institute of Health Rapid Acceleration of Diagnostics Program.

