

# Forward-Looking Statements & Pro-Forma 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 proforma financial measures. Management uses these pro-forma 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 other period results and are useful to investors and financial analysts in assessing the Company's operating performance. The pro-forma 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 pro-forma measures to their most directly comparable GAAP financial measures set forth herein and in the earnings press release issued earlier today.



#### **Progress on Assay Redevelopment Roadmap**

Q2 2022

- CEO transition on April 25th
- Business strategic review, starting in June



Q3 2022

- Strategic review completed in July, business realignment and restructure in August
- Re-development team and charter established
- Shelf-life of our products phase 1



Q4 2022

- Stability and variance improve ments for common assay components
- Identify primary wave of improvements transitioned into manufacturing



- Manufacturing implementation of primary wave
- Raw material specifications, stability and qualification
- Automated WIs & Documentation



- Process automation
- Process scaling
- Optimized workflow
- Product and process harmonization
- Shelf-life
   Enhancements

28.3%

Pro-forma Gross Margin\*

Revenue

\$23.5M

\$26.6M

35.0%

\$24M-\$26M

\*Proforma reconciliation included on Slide 6

**Quanterix** 

... Progress according to plan

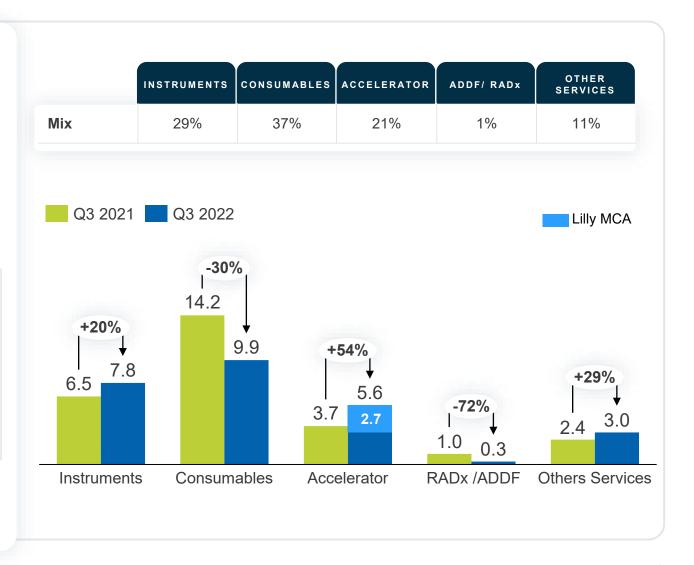
## Q3 2022 Revenue results are aligned with company expectations

Reiterate full year 2022 revenue to be flat compared to 2021 \$105M (excluding one time '21 NIH grant funding)

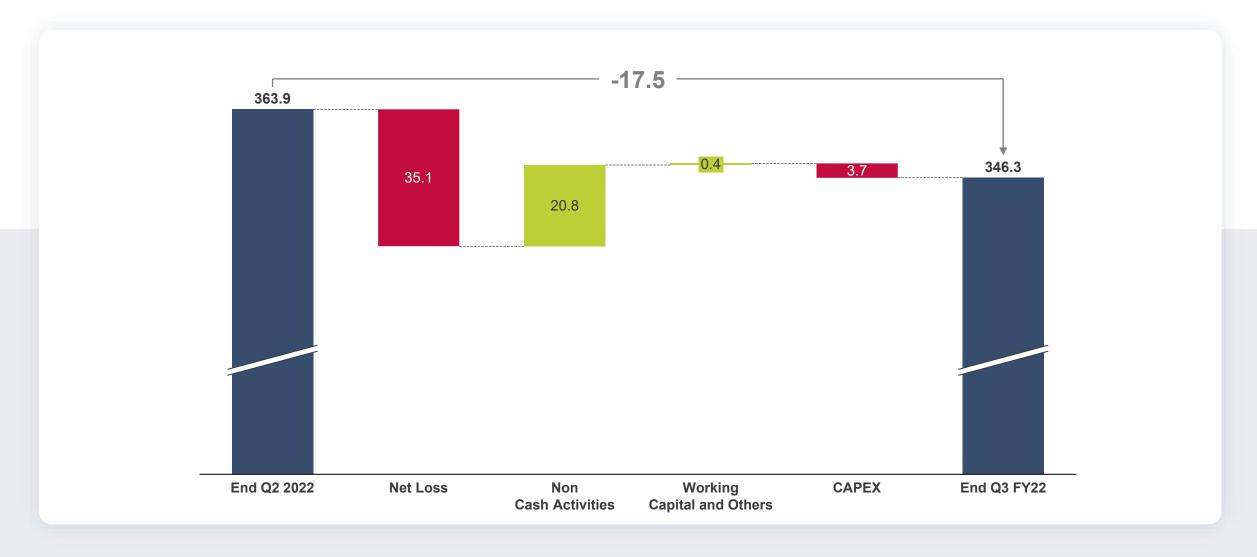
	Q3 GA	AAP	Q3 Proforma					
	2021	2022	2021	2022				
Revenue	27.7 (26.7 excl. RADx)	26.6	27.7 (26.7 excl. RADx)	26.6				
GM \$	15.2	10.9	13.8	9.3				
GM %	55.1%	41.1%	49.8%(1)	35.0%(1)				
Operating Expense	30.5	26.6	29.0(1)	25.0(1)(2)				
Operating Loss	-15.2	-15.7	-15.2	-15.7				
Operating Loss as % of Revenue	-55.0%	-58.7%	-55.0%	-58.7%				

<sup>(1)</sup> Proforma reconciliation included on Slide 6.

<sup>(2)</sup> Excludes restructuring and impairment expenses incurred in Q3 2022 in connection with our restructuring plan. See Supplemental Schedule of Operating Expenses on Slide 7 for reconciliation.



#### Q3 2022 Cash & Cash Burn





#### Reconciliation of GAAP to Pro Forma

# Quanterix Corporation Reconciliation of GAAP Financial Measures to Non-GAAP Financial Measures (Unaudited and in thousands, except percentages)

		Three Months Ended September 30,				Three Months Ended June 30,		Nine Months Ended September 30,			
		2022		2021		2022		2022		2021	
GAAP gross profit	\$	10,944	\$	15,244	\$	8,711	\$	34,214	\$	45,467	
Distributions costs (1)		(1,617)		(1,462)		(2,065)		(5,546)		(5,024)	
Non-GAAP pro forma gross profit	\$	9,327	\$	13,782	\$	6,646	\$	28,668	\$	40,443	
Gross margin		41.1%		55.1%		37.1%		42.9%		56.6%	
Non-GAAP pro forma gross margin		35.0%		49.8%		28.3%		36.0%		50.4%	
GAAP total operating expenses  Distributions costs (1)  Non-GAAP pro forma total operating costs	\$ 	47,547 (1,617) 45,930	\$ 	30,477 (1,462) 29,015	\$ 	33,670 (2,065) 31,605	\$ 	113,963 (5,546) 108,417	\$ 	84,157 (5,024) 79,133	
GAAP loss from operations	<u>\$</u>	(36,603)	<u>\$</u>	(15,233)	<u>\$</u>	(24,959)	<u>\$</u>	(79,749)	\$	(38,690)	
Non-GAAP pro forma loss from operations	<u>\$</u>	(36,603)	\$	(15,233)	<u>\$</u>	(24,959)	<u>\$</u>	(79,749)	<u>\$</u>	(38,690)	

<sup>(1)</sup> Distribution costs, which include freight and other activities costs associated with product shipments, net of charges passed on to the customer, are captured within operating expenses in our consolidated statements of operations. During the three and nine months ended September 30, 2022, we incurred \$1.6 million and \$5.5 million, respectively, of distribution costs recorded within operating expenses. During the three and nine months ended September 30, 2021, we incurred \$1.5 million and \$5.0 million, respectively, of distribution costs recorded within operating expenses. During the three months ended June 30, 2022, we incurred \$2.1 million of distribution costs recorded within operating expenses.



# Supplemental Schedule of Operating Expenses

Quanterix Corporation
Supplemental Schedule of Operating Expenses
(Unaudited and in thousands)

		Three Months End	ded Se	eptember 30,		Nine Months End	ded September 30,		
		2022		2021		2022		2021	
Operating expenses:									
Research and development	\$	6,631	\$	6,807	\$	20,290	\$	20,244	
Selling, general and administrative		19,966		23,670		72,723		63,913	
Subtotal	\$	26,597	\$	30,477	\$	93,013	\$	84,157	
Other lease costs		609				609			
Restructuring		3,426		_		3,426		_	
Goodwill impairment		8,220		_		8,220		_	
Impairment expense		8,695				8,695			
Subtotal	\$	20,950	\$	<u> </u>	\$	20,950	\$	<u> </u>	
Total operating expenses	<u>\$</u>	47,547	<u>\$</u>	30,477	<u>\$</u>	113,963	\$	84,157	

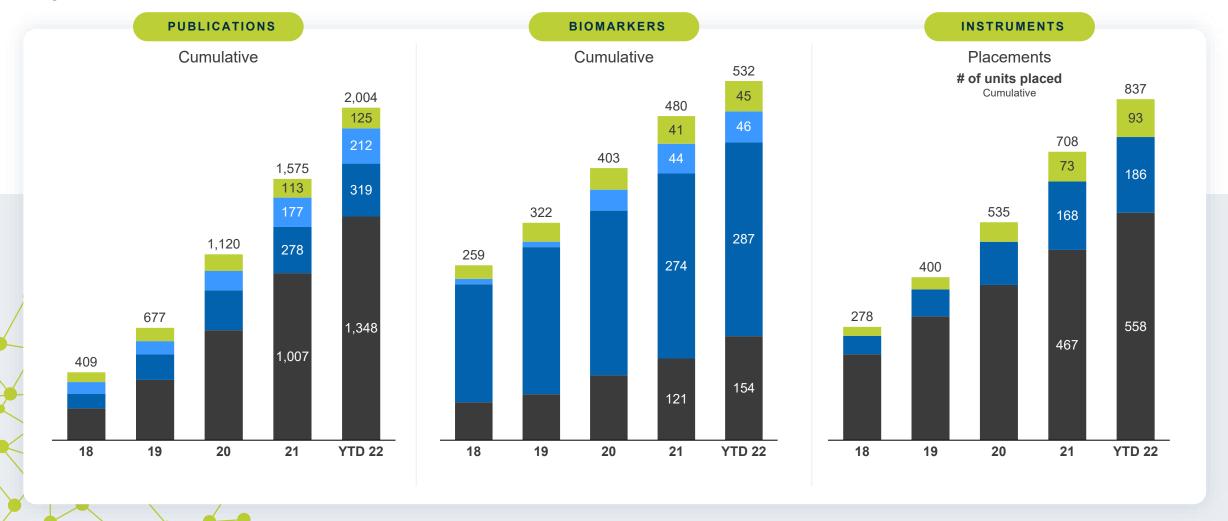
	Three Months Ended September 30,					Nine Months Ended September 30,				
	2022			2021		2022	2021			
Operating expenses:										
Research and development	\$	6,631	\$	6,807	\$	20,290	\$	20,244		
Selling, general and administrative		19,966		23,670		72,723		63,913		
Subtotal	\$	26,597	\$	30,477	\$	93,013	\$	84,157		
Distributions costs (1)		(1,617)		(1,462)		(5,546)		(5,024)		
Subtotal proforma OpEx	\$	24,980	\$	29,015	\$	87,467	\$	79,133		
Other lease costs		609		_		609		_		
Restructuring		3,426		_		3,426				
Goodwill impairment		8,220				8,220				
Impairment expense		8,695				8,695				
Subtotal	\$	20,950	\$		\$	20,950	\$			
Non-GAAP pro forma total operating costs	<u>\$</u>	45,930	\$	29,015	\$	108,417	\$	79,133		



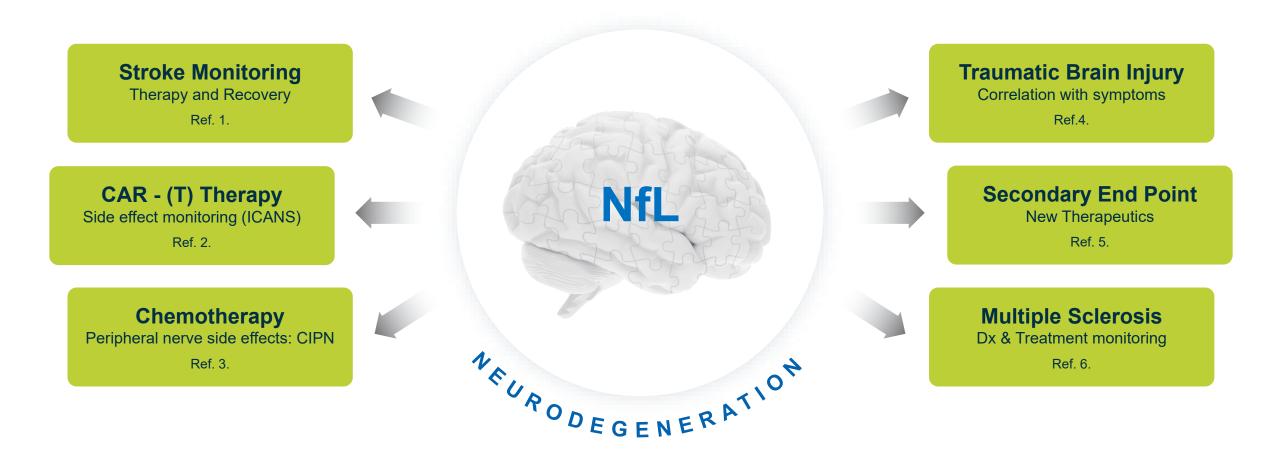
# Scientific Validation Driving Adoption

2022 Advances





#### New Simoa NfL Publications: Key Biomarker to Assess Neural Damage



- 1. Wang et.al., Front. Neurol., 25 October, 2022; 2. Butt et.al., JAMA Oncology, 1 September, 2022; 3. Brugess et.al., Nature, 16 September, 2022
- 4. Lange et.al., J. Neurotrama, 18 October, 2022; 5. Imamura et.al., J,eClinn. 17 October, 2022; 6. Williams et.al., Mult. Scler. 28 October, 2022

## Alzheimer's: Blood Biomarkers Increasingly Utilized in Clinical Trials

#### **Accelerants for efficiency & effectiveness:**

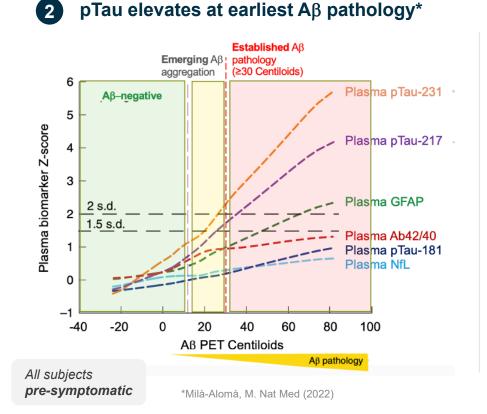
- 1 Pre-screening/enrollment of larger, more diverse populations with decrease screen fails 3 Monitoring of drug effects (pharmacodynamic)
- 2 Improved targeting along the disease continuum; entry criteria for preventative AD trials
- 1 pTau reduces PET screen fails\*

Proportion of pts Aβ and Tau PET positive using plasma pTau-181 as a pre-screener

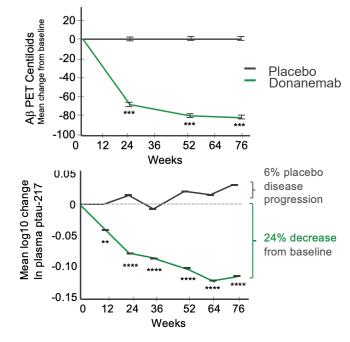




■ PET positive ■ PET negative



#### 3 pTau tracks w/ Aβ removal



\*TRAILBLAZER-2 data, presented at AAIC 2021