

Corporate Presentation

August 2024

Quanterix™
Discovery Fueled by Ultra-Sensitivity



Forward-Looking Statements & Non-GAAP Financial Measures

This presentation contains forward-looking statements within the meaning of the U.S. 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's 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's actual results to differ from those expressed or implied in the forward-looking statements in this presentation are discussed in Quanterix's 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 Quanterix's financial statements presented on a GAAP basis, Quanterix has provided certain non-GAAP financial measures. Management uses these non-GAAP measures to evaluate our operating performance in a manner that allows for meaningful period-to-period comparison and analysis of trends in our business and our competitors. Management believes that presentation of these non-GAAP measures provides useful information to investors in assessing our operating performance within our industry and in order to allow comparability to the presentation of other companies in our industry. 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 herein and in the associated earnings press release.

Unless otherwise specified, all financial information contained herein is provided as of June 30, 2024, except that all guidance is as of August 8, 2024. This document is not a reaffirmation of guidance.



Quanterix[®]

Our Mission

Create the tools enabling discovery and better health

Strong Research Business

Business model redefined to enable innovation, expected to grow double-digits¹

Differentiated Technology

The power of ultra-sensitive biomarker detection with Simoa[®]

Diagnostics Opportunity

Built to lead the Neuro Decade



Quanterix®

Strong Research Business

Business model redefined to enable innovation, expected to **grow double-digits**¹

Quanterix Today



Unmatched Technology



~1,000 Instruments installed



Early penetration in multi-billion diagnostics



CLIA lab running validated neuro LDTs

2024 Guidance¹

\$136M

Revenue

53%

Non-GAAP
Gross Margin²

\$30M

Cash burn

Q2'24 Ending

\$300M

Cash Balance³

560+

Biomarkers

2,900+

Publications

 Nasdaq **QTRX**

1. Full year 2024 Guidance at the mid-point for Revenue and Non-GAAP gross margin. The Company has guided cash burn of \$30M for full year 2024.
2. The Company has guided full year 2024 GAAP gross margin percentage to be 57-61%, and Non-GAAP gross margin percentage to be 51-55%. GAAP gross margin does not include shipping and handling costs, which include freight and other activities costs associated with product shipments. Non-GAAP gross margin includes these shipping and handling costs.
3. Includes Cash, Cash Equivalents, Marketable Securities, and Restricted Cash Balances.

Integrated Business Model, Buffer Against Macro Headwinds

With ~80% of recurring revenue mix in Q2'24¹

DISCOVERY TO DIAGNOSTICS

Discovery

Instruments

- Feature both bead assay and planar array technology
- Used in 2,900+ publications



Consumables & Reagents

- A broad selection of assays developed for a wide range of therapeutic areas

Clinical Trials

Accelerator Services

- Contract research services through Simoa Accelerator Laboratory
- Sample Testing Services
- Custom Assay Development

Diagnostics

CLIA Diagnostic Lab

- Enabling the future of Therapeutics for Neurodegenerative Diseases
- p-Tau 217 LDT
- p-Tau 181 LDT
- Nf-L LDT

Accelerator Lab

A unique and durable model, with a full suite of services for exploratory research, clinical trials, and diagnostics¹



Expertise

150+ years of
assay development
experience²

Supporting biomarker
research and clinical trials



Product Incubator

2300+ projects
and **480+**
customers globally

Insights into key biomarkers,
leading to assay development



Commercial Engine

37%
revenue growth³

Valuable offering, offsetting
weak instrument environment

Efficient operating framework in place

Improvement of **50%**
in 6 quarters¹

Testing Scale

**>0.75M tests
per year**

with capacity to
increase to 3x

**Accelerator
Lab**

Improvement of **300%**
in 6 quarters¹

Manufacturability

**>4M tests
per year**

with capacity to
increase to 3x

**Assay
Manufacturing**

Increase of **80%**
in 6 quarters¹

Studies

**>45 ongoing
studies**

supported in
Neurology

**Research
partnerships**

Strong foundation for a ramping innovation rate in 2024



Prior 2 years

2023

2024 and beyond

Launched **0**
new assays

One LDT Test

Launched
5 new assays

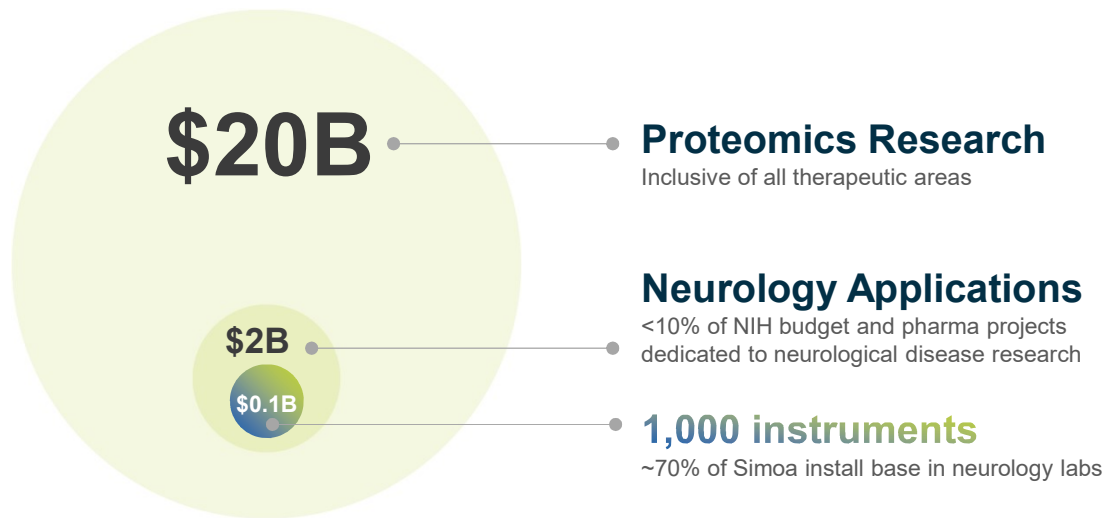
One LDT Test

≥20 assays a year
Develop platform
(Plexity & automation)
IVD launch

Harmonized processes driving rapid menu expansion

Assay development times reduced from >18 months to <6 months

Simoa[®] - in every research lab



Opportunities to unlock TAM

Therapeutic Area Expansion

Growth into new applications and biomarker assays to capture research market outside of neurology

Increased Sensitivity

Continued innovation in driving deeper levels of insight by enabling further advancement in ultra-sensitive quantitative proteomics

Increased Access

Investment in future solutions to democratize Simoa platform (footprint, price, workflow)



100,000+

Research Labs:

- Expanded therapeutic areas
- International
- Government institutions
- Private institutions



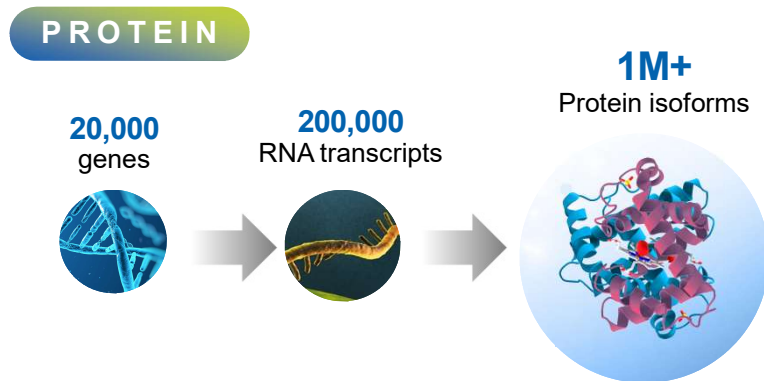
Quanterix[®]

Differentiated Technology

The power of ultra-sensitive biomarker detection with Simoa[®]

Proteome is dynamic and closest to clinical actionability

Simoa[®] sensitivity allows digital detection of proteins and their isoforms



Simoa[®] sensitivity enables...

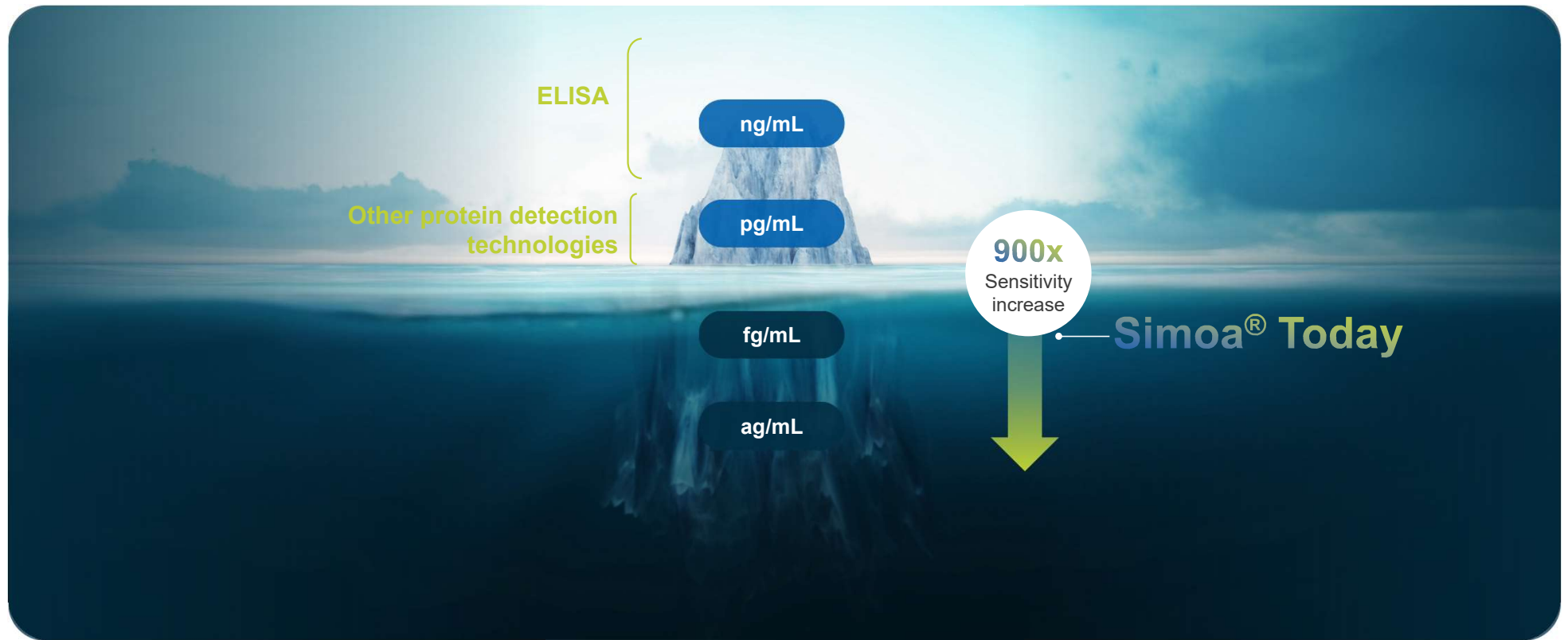
Less invasive, smaller sample

New biomarkers & protein isoforms

Multiplexing for disease specificity

Single molecule to unravel heterogeneity

Simoa[®] detects biomarkers unmeasurable by other technology

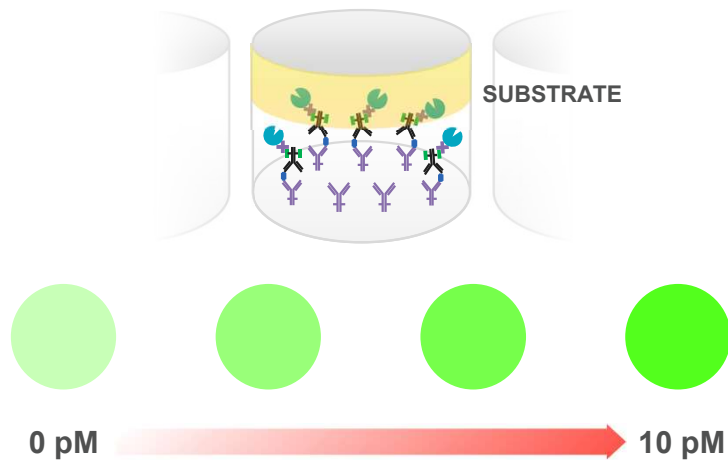


Surpassing detection limits to deliver exquisite sensitivity

Analog

Traditional ELISA assay

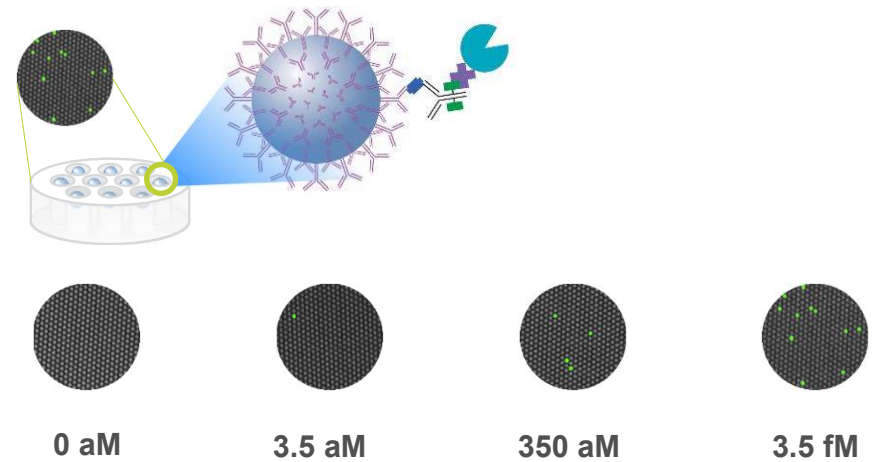
Millions of molecules needed to reach detection limit



Simoa[®]

Digital

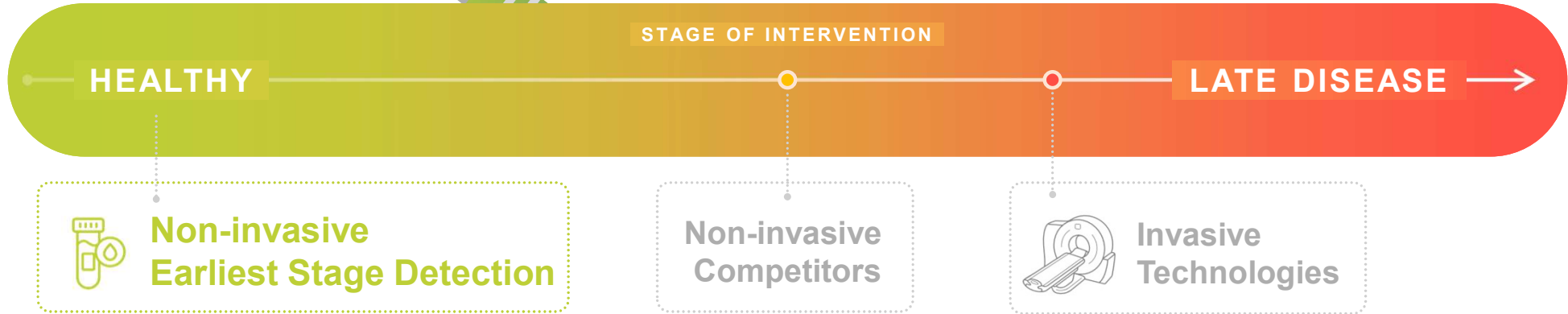
Single molecule needed to reach detection limit





Catching disease early enables health-care...

...versus today's sick-care management





Quanterix[®]

Diagnostics Opportunity

Built to lead the Neuro Decade

The decade of Neuro





Source:

1. 2023 Alzheimer's Disease Facts and Figures from Alzheimer's Association® / Alzheimer Disease International (<https://www.alzint.org>)

2. <https://clinicaltrials.gov/> and company reports

Multiple ongoing studies in Neurology

- ✓ AD
- ✓ ALS
- ✓ MS
- ✓ Parkinson
- ✓ TBI
- ✓ Neuropsychology



**With
Alzheimer's Disease
as the core of our
Diagnostics focus**

Alzheimer's Diagnostics Opportunity



>55M globally
living with Alzheimer's

By 2050 projected to
rise to 139M

TD TD Cowen
a division of TD Securities

\$3B

LEERINK PARTNERS

\$9B

Goldman
Sachs

\$10B

A significant TAM¹ to
unlock

Drivers

Therapeutic Uptake

Reimbursement

Regulatory Approval



Mortality in the US more than breast and prostate cancer, combined.



¹TAM estimated for blood-based biomarker testing in Alzheimer's diagnostics.

Early detection of Alzheimer's pathology requires greater sensitivity

7M with AD in US

Addressable by

LIMIT OF DETECTION OF OTHER IMMUNOASSAYS

>140M over the age of 45 in U.S.

25M w/ early amyloid prevalence^(b)

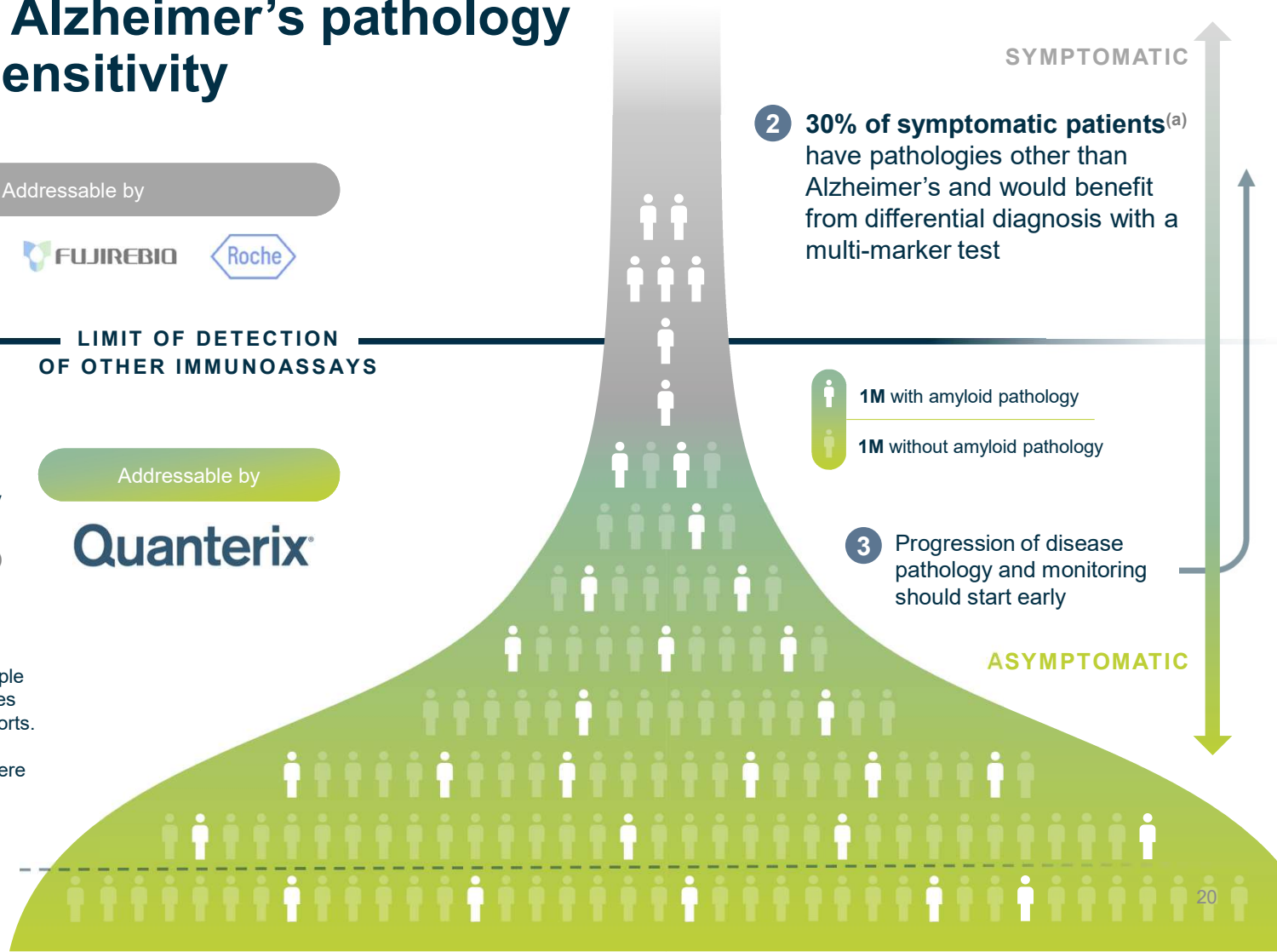
Addressable by

1 Quanterix has a single femtogram/mL limit of detection and has yet to encounter a single sample unreadable below this level in over 2,000 samples tested, including the BioHermes and VUMC cohorts. In similar patient cohorts, other technology platforms have shown up to ~30% of samples were unreadable below the platform detection limit.

FUTURE CLINICAL CUT OFF

Quanterix
Discovery Fueled by Ultra-Sensitivity

Source:
a. WHO
b. JAMA Neurology



2 30% of symptomatic patients^(a) have pathologies other than Alzheimer's and would benefit from differential diagnosis with a multi-marker test

1M with amyloid pathology
1M without amyloid pathology

3 Progression of disease pathology and monitoring should start early

Simoa®: A Result for Every Patient

SIMOA

1000 patients



ALL
patients
receive
a result¹

Mass Spectrometry

1000 patients



280
patients
with NO
result²



Chemiluminescence

1000 patients



390
patients
with NO
result³



While most p-Tau 217 tests can achieve high accuracy (e.g., AUC), the best clinical utility is when all patients receive a result.

1. With its single femtogram/mL limit of detection, SIMOA has provided p-Tau 217 results in all patient samples tested, including over 2,000 samples in BioHermes and VUMC cohorts.

"Clinical validation of the LucentAD p-Tau 217 as a lab developed test (LDT) for clinical use." Presented at AD/PD, March 2024.

"Multi-marker approach to reducing the intermediate range of a high accuracy 2-cutoff plasma p-Tau 217 test for amyloid detection." Presented at AAIC, July 2024.

<https://www.quanterix.com/whitepapers-appnotes/>

2. June 2024: FNIH head-to-head study: *"Of the 392 samples included, 110 samples had p-tau217 values that were below the level of detection for the C2N PrecivityAD2 assay..."*

"Notably, both the C2N PrecivityAD2 and Fujirebio Lumipulse p-tau217 assays had limited sensitivity, resulting in imputed values for samples with low values..."

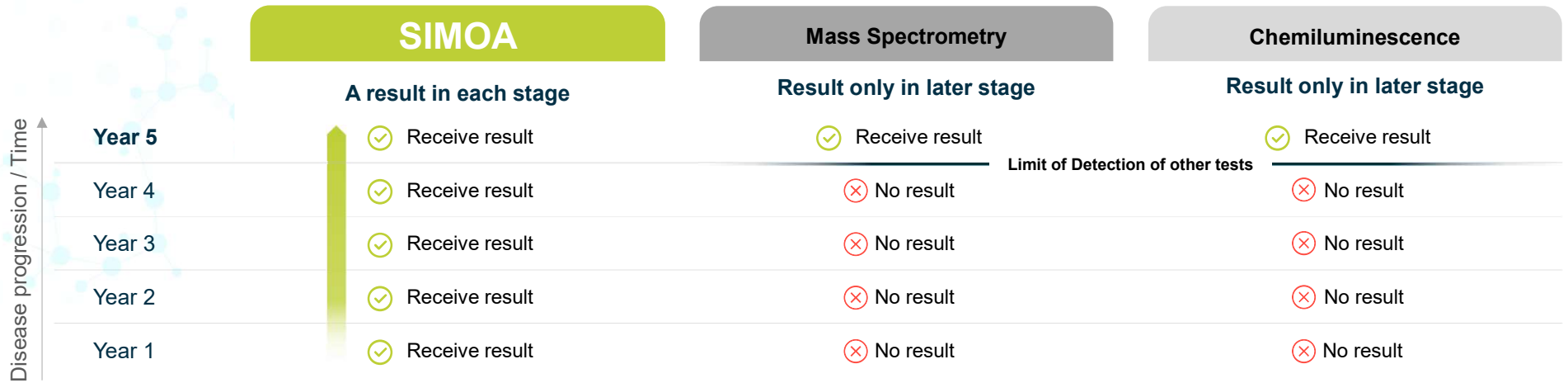
SIMOA read every patient sample in the same cohort.

[FNIH head-to-head study](#)

3. August 2022: Roche in Alzheimers Dement: *"The main limitation of the present study was that measured plasma P-tau217 N-terminal and mid-domain concentrations were in many cases below the lower level of detection. For P-tau217 N-terminal, 39% had to be assigned the calculation minimum level..."*

[Alzheimers Dement](#)

Simoa[®]: Best in class for monitoring early AD




Family history or Memory concerns → Visit neurologist

Get a result
 Quantifiable Tau level result each year to track progression and potentially earlier treatment

Get no result

Physician may need to use other tests to follow up e.g., PET, CSF or SIMOA.

Quanterix best positioned for Alzheimer's detection today

 **Quanterix**
Enabled DX lab

 **Non-Quanterix**
DX lab

Simoa[®]

Mass Spectrometry

Chemiluminescence

**Quantitative range
in clinical cohorts**

High

High

Limited

**Detect p-Tau 217
(Meet NIA-AA guidelines)**

Yes

Yes

No

Cost of test

Low

<\$400/test

High

>\$1200/test

Low

<\$400/test

Scalable platform

Yes

No

Yes

**Multiplex capabilities
(for differentiated DX)**

Yes

Limited

No

**Simoa[®] technology
uniquely suited for
building global
Diagnostics testing
infrastructure**

Multi-marker: the Next Phase of AD Testing

Today: Single Marker p-Tau 217



**p-Tau 217 immunoassays
launched or in development**

Tomorrow: Multi-Marker Test for AD

Quanterix[®]
**p-Tau 217, A β 40/42,
NfL, GFAP**

A better test, with an algorithm, and
potential for differential diagnosis

✓ **Accurate**
+90% Sensitivity
+90% Specificity
+90% Accuracy

✓ **Actionable**
10.5% Indeterminate vs.
>30% Indeterminate for
stand-alone p-Tau 217

Expect to launch Multi-Marker Test by end of 2024

Alzheimer's Diagnostic Investment and Execution

Allocating capital

Over \$20M¹ of capital
allocated for next 2 years to
**advance AD
Diagnostics**

to execute commercial adoption

Reimbursement:

CPT codes (Pricing expected later in 2024)

FDA:

Breakthrough designation received for 3 tests

Updates on Key Studies:

CANTATE: P1 Complete, P2 in progress, P3 Staged

BioHermes: P1 Complete

Multi-marker data presented at AAIC in July 2024

¹\$10-15M of this investment expected in 2024.

**Diversified portfolio
geared to beat industry
headwinds with recent
transformation and
multiple growth drivers**

Investment Thesis

- ✓ Completed transformation provides **leverage for growth** and expanding gross margins
- ✓ **Strong balance sheet**, with no near-term financing needs
- ✓ A **diverse, resilient business model**, indexed to recurring revenue and growing double digits
- ✓ **Differentiated technology** moat driving Research and new exciting opportunities in Diagnostics



Appendix

Executive Team



Chief Executive Officer
Masoud Toloue
Revvity, GenoHub, Bioo Sc.



Chief Financial Officer
Vandana Sriram
Azenta, GE



Chief People Officer
Erica Bell
BICO, Revvity, GE



SVP Corp Dev. & Strategy
Naren Bhat
Revvity, GE, Philips



SVP & General Counsel
Laurie Churchill
LeMaitre Vascular, Avid Tech,
Ropes & Gray



Chief Commercial Officer
Darrin Crisitello
Mission Bio, Natera, Color
Health



Chief Information Officer
Alexandra Phillips
Hologic, Revvity, Boston
Scientific, Tyco



Chief Operating Officer
Mike Miller
ProterixBio, BioScale, Axsun
Tech, Physical Sciences Inc.



**Chief Science and
Collaboration Officer**
Mark Roskey
Revvity, Caliper Life Sc.

Products and Services

Simoa® Ultra-sensitive Instruments



HD-X

Bead assay technology

Multiplexing up to 4-plex

Floor-standing integrated automated assay prep and detection, *sample in, answer out*



SR-X

Bead assay technology

Multiplexing up to 4-plex

Benchtop form factor semi-automated offline assay prep



SP-X

Planar array technology

Multiplexing up to 10-plex

Benchtop form factor semi-automated offline assay prep

Assay Kits



Bead-based & Planar array

Broad selection of catalog and custom assays developed for neurology, oncology, cardiology, infectious disease and immunology research

Capable of assay customization with homebrew kits

Single-plex and multiplex formats

Services



CRO

Contract research services through Simoa Accelerator Laboratory

Sample testing services

Custom assay development

Custom reagent production and kitting

Diagnostics

CLIA – certified lab

Menu of Laboratory Tests available

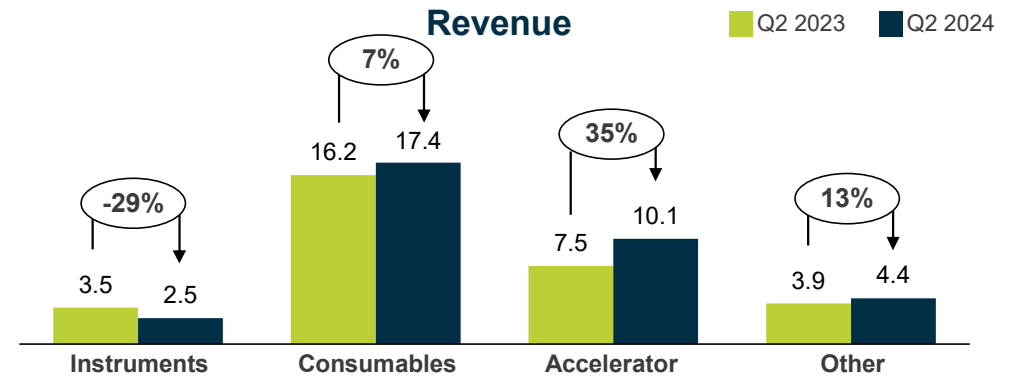
Direct to Patient and Provider portal

Q2 2024 Results vs PYQ2

(in millions)

	Q2 GAAP		Q2 Non-GAAP	
	2023	2024	2023	2024
Revenue	31.0	34.4	31.0	34.4
Gross Margin \$	19.1	20.1	17.5	18.0
Gross Margin %	61.7%	58.3%	56.4%	52.3%
Operating Expense	28.7	33.2	27.1	31.1
Operating Loss	-9.6	-13.1	-9.6	-13.1
Cash Usage	-0.1	-5.1	-0.1	-5.1

	INSTRUMENTS	CONSUMABLES	ACCELERATOR	OTHER
Revenue Mix	7%	51%	29%	13%

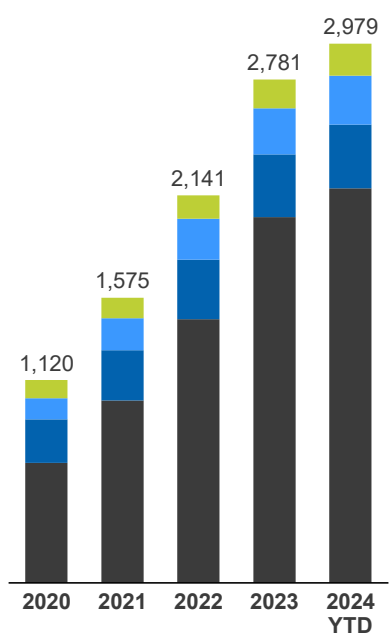


Scientific Validation Driving Adoption



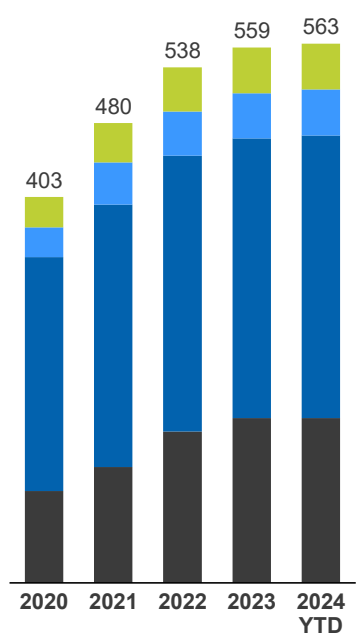
PUBLICATIONS

Cumulative



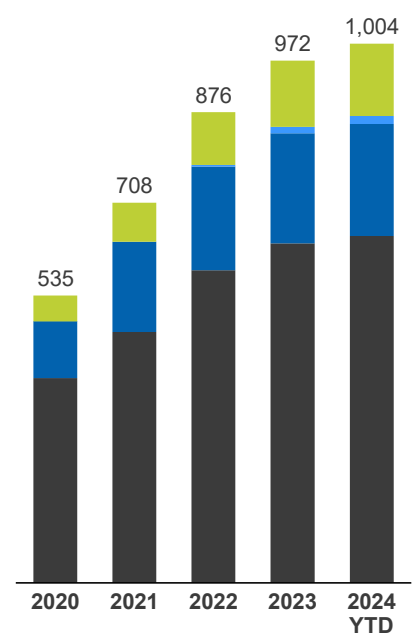
BIOMARKERS

Cumulative



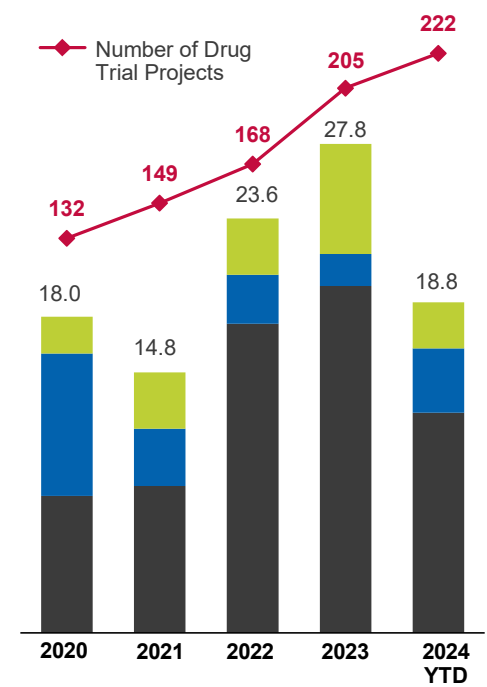
INSTRUMENTS

Placements
of units placed
Cumulative



ACCELERATOR

Projects & Revenue



Reconciliation of GAAP to Non-GAAP Financial Measures

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

	Three Months Ended June 30,		Six Months Ended June 30,	
	2024	2023	2024	2023
GAAP gross profit	\$ 20,058	19,138	\$ 39,684	\$ 36,064
Shipping and handling costs (1)	(2,075)	(1,623)	(4,217)	(3,451)
Non-GAAP gross profit	\$ 17,983	17,515	\$ 35,467	\$ 32,613
GAAP revenue	\$ 34,381	31,029	\$ 66,447	\$ 59,485
GAAP gross margin (gross profit as % of revenue)	58.3%	61.7%	59.7%	60.6%
Non-GAAP gross margin (non-GAAP gross profit as % of revenue)	52.3%	56.4%	53.4%	54.8%
GAAP total operating expenses	\$ 33,166	28,699	\$ 66,758	\$ 55,045
Shipping and handling costs (1)	(2,075)	(1,623)	(4,217)	(3,451)
Non-GAAP total operating expenses	\$ 31,091	27,076	\$ 62,541	\$ 51,594
GAAP loss from operations	\$ (13,108)	(9,561)	\$ (27,074)	\$ (18,981)
Non-GAAP loss from operations	\$ (13,108)	(9,561)	\$ (27,074)	\$ (18,981)

(1) Shipping and handling costs, which include freight and other activities costs associated with product shipments, are captured within operating expenses in our consolidated statements of operations.