

J.P. Morgan Healthcare Conference 2024

Masoud Toloue, CEO

January 10, 2024

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'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.

Forward-looking statements in this presentation include certain preliminary, unaudited estimated financial information for the fourth quarter and year ended December 31, 2023. This financial information is unaudited and preliminary and does not present all information necessary for an understanding of Quanterix's financial condition as of December 31, 2023 and its results of operations for the quarter and year ended December 31, 2023. This financial information is subject to completion of normal quarter and year-end close procedures. These procedures and the audit of Quanterix's financial statements for the year ended December 31, 2023 are ongoing and could result in changes to this financial information.

To supplement Quanterix's financial information 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 as a factor in assessing progress against our corporate transformation. 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 corresponding GAAP financial measures.

Create the tools enabling discovery and better health



Unmatched Technology



>950 Instruments installed



Early penetration in multi-billion diagnostics



CLIA lab running validated neuro LDTs

\$121M

Revenue¹

\$18M

Cash burn¹

High **50%+'s**

GAAP
Gross Margin²

Appr. **50%**

Non-GAAP
Gross Margin²

550+

Biomarkers

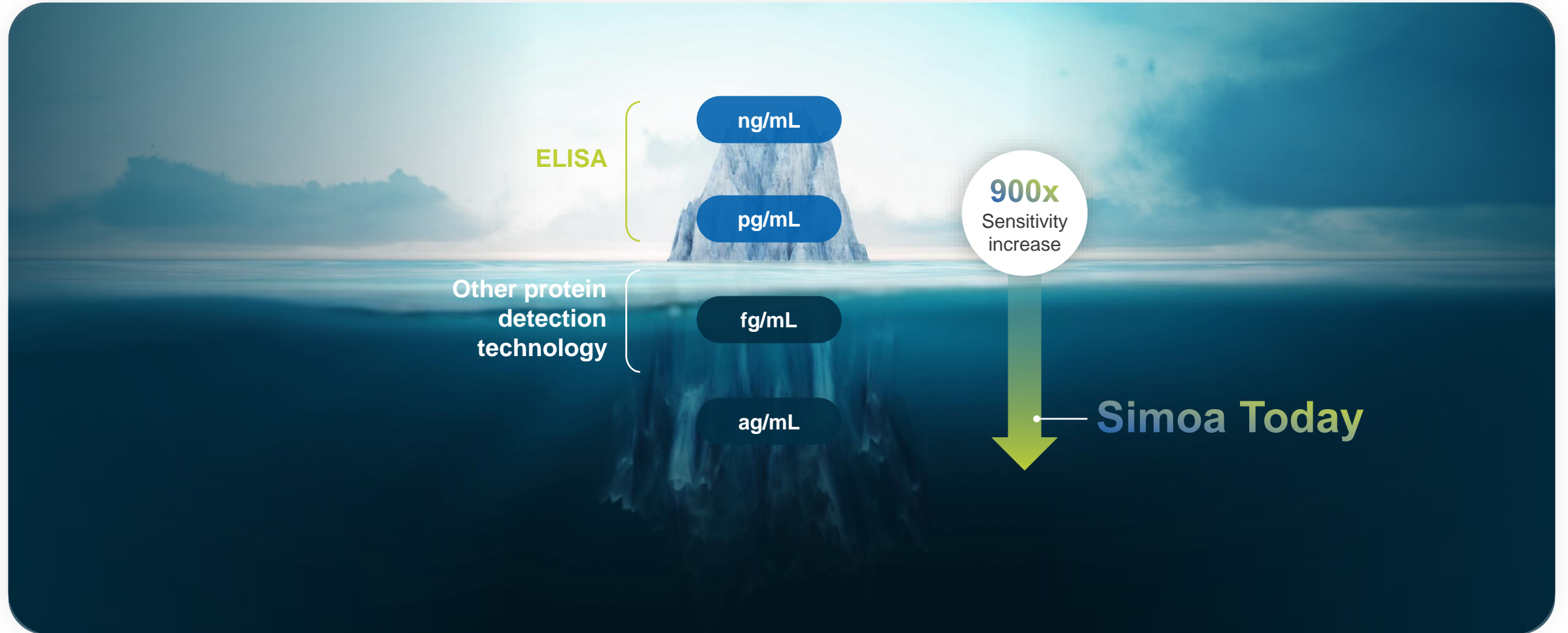
2,500+

Publications

 Nasdaq : QTRX

1. Preliminary unaudited full year results for 2023.
2. The Company is guiding full year 2023 GAAP gross margin percentage to be in the high 50's, and non-GAAP gross margin percentage to be approximately 50%. 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.

Simoa detects biomarkers unmeasurable by other technology

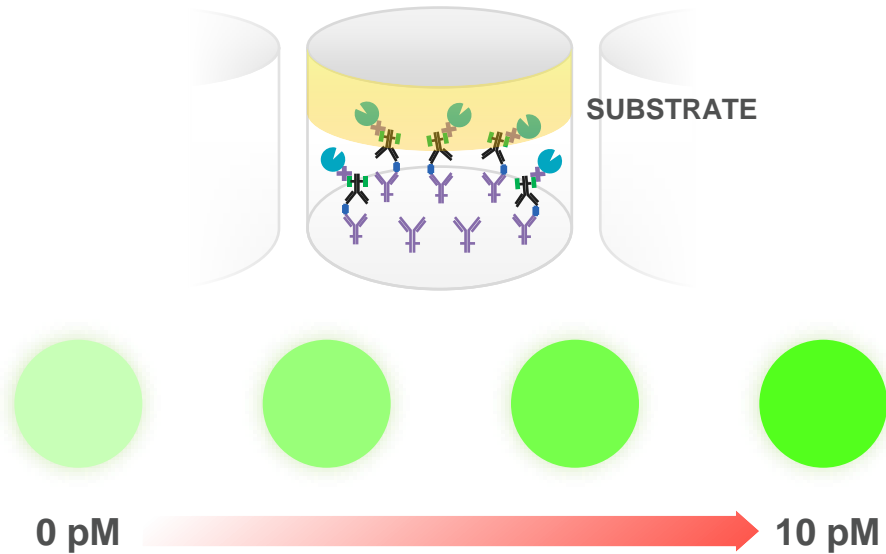


Surpassing detection limits to deliver exquisite sensitivity

Analog

Traditional ELISA assay

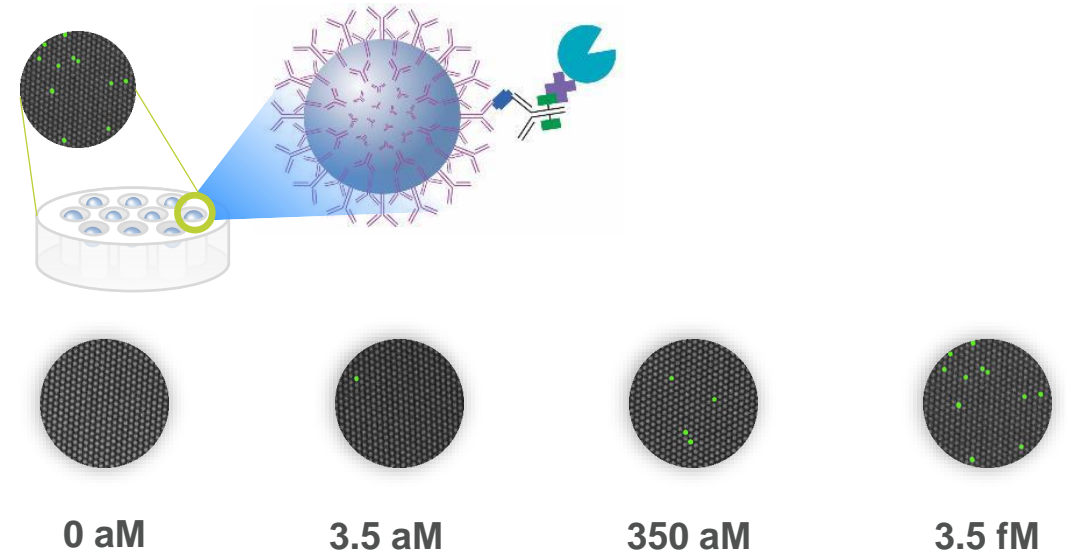
Millions of molecules needed to reach detection limit



Simoa[®]

Digital

Single molecule needed to reach detection limit





Catching disease early enables health-care...

...versus today's sick-care management

STAGE OF INTERVENTION

HEALTHY

LATE DISEASE

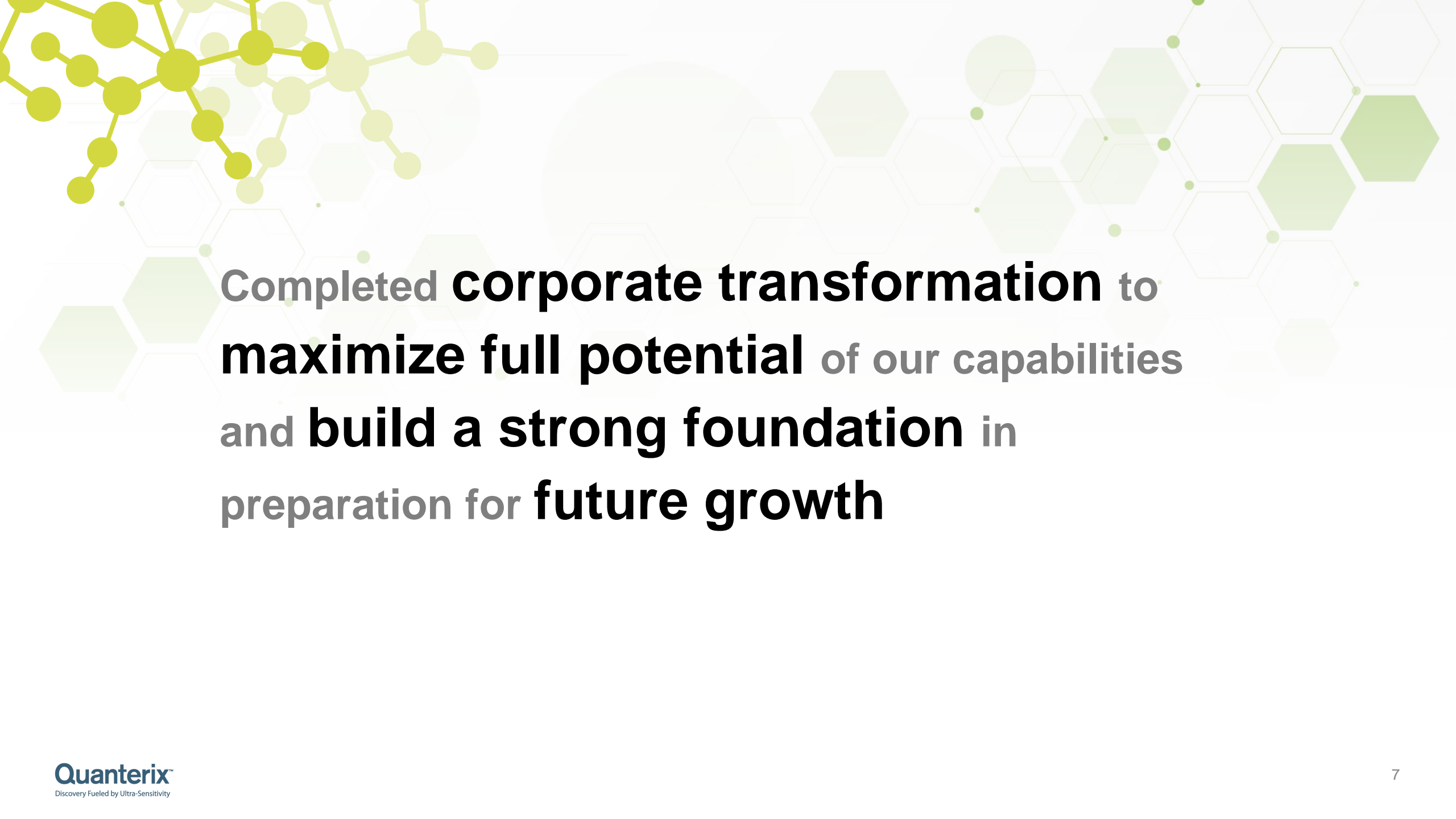


**Non-invasive
Earliest stage detection**

Non-invasive
Competitors

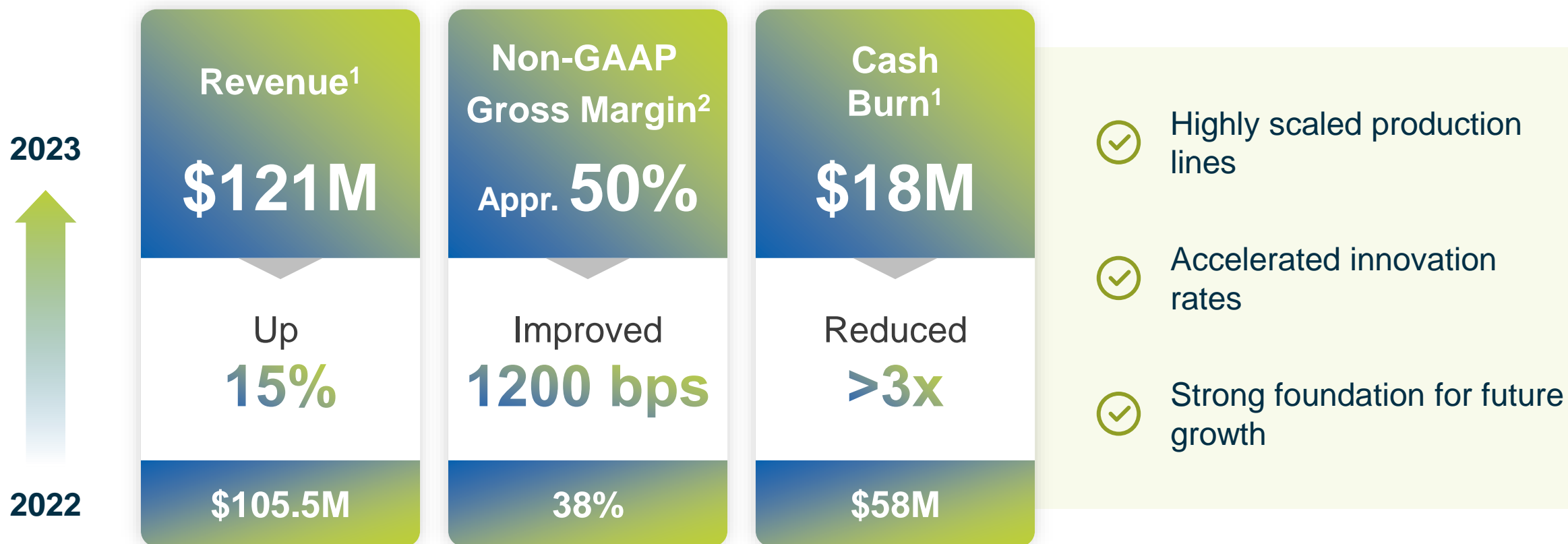


Invasive
Technologies



Completed **corporate transformation** to
maximize full potential of our capabilities
and **build a strong foundation** in
preparation for **future growth**

Consistent execution while driving broad transformation



1. Preliminary unaudited full year results for 2023.

2. The Company is guiding full year 2023 GAAP gross margin percentage to be in the high 50's, and non-GAAP gross margin percentage to be approximately 50%. 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.

Delivering an efficient operating framework

Improvement of **50%**
in 6 quarters

Testing Scale

**>0.75M¹ tests
per year**

with capacity to
increase to 3x

**Accelerator
Lab**

Improvement of **300%**
in 6 quarters

Manufacturability

**>4M¹ tests
per year**

with capacity to
increase to 3x

**Assay
Manufacturing**

Increase of **80%**
in 6 quarters

Studies

**>45 ongoing
studies**

supported in
Neurology

**Research
partnerships**

As the foundation for a ramping innovation rate in 2024



Prior 2 years

Launched **0**
new assays

One LDT Test

2023

Launched **5**
new assays

One LDT Test

2024 and beyond

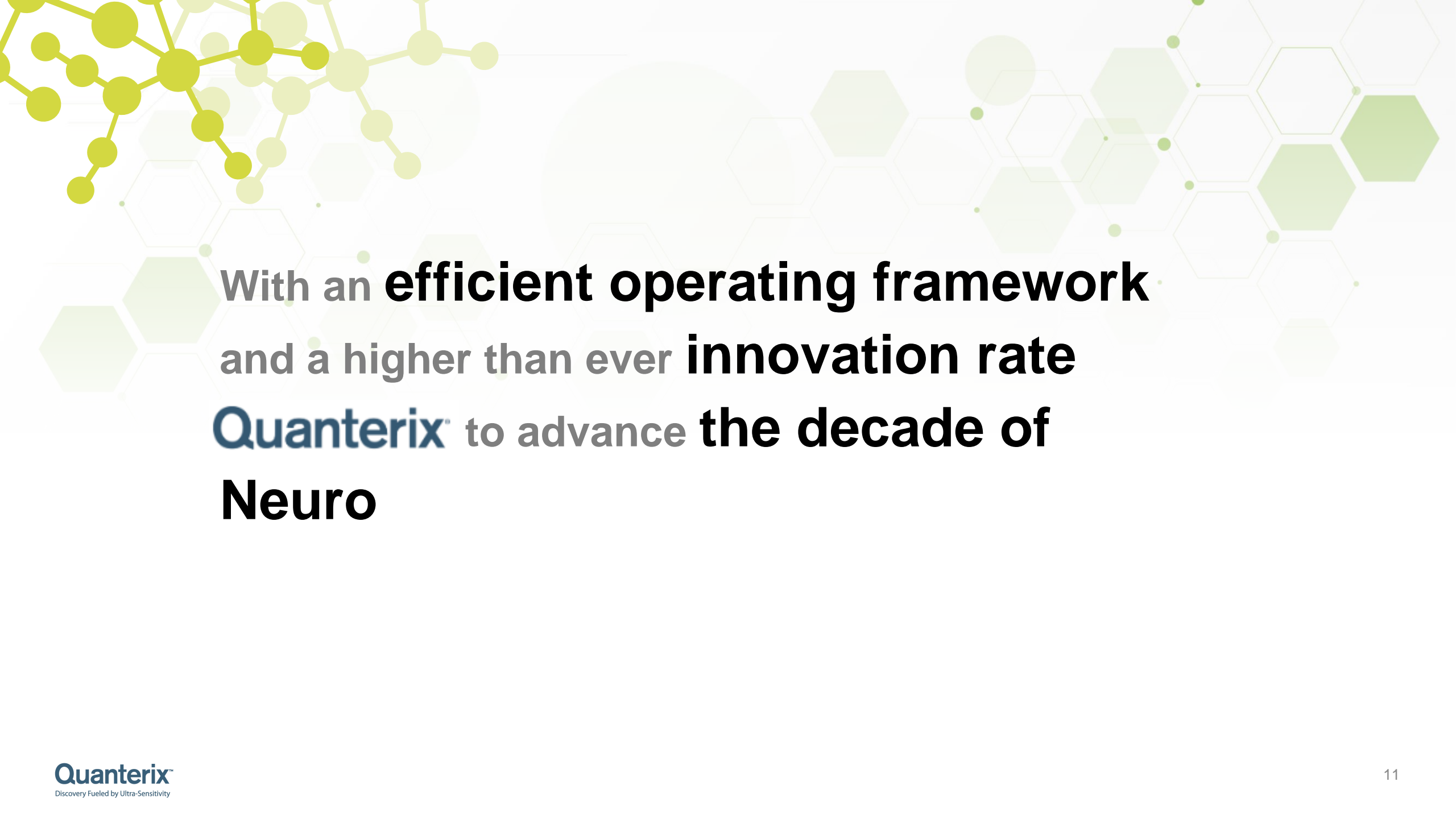
>20 assays a year

Develop platform
(Plexity & automation)

IVD launch

Harmonized processes driving rapid menu expansion

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



With an **efficient operating framework**
and a higher than ever **innovation rate**
Quanterix to advance **the decade of**
Neuro

The decade of Neuro

Increasing Demographic Need

Projected **139M AD cases**¹ by 2050

Extensively validated blood biomarkers

Supporting FDA **accelerated approval** (e.g. NfL)

Approved therapies

First drugs approved in 2023
EISAI / Biogen



Emerging pipeline

>20 large biopharma
running >120 CNS programs²

Accelerator Lab indicators

47 neuro trials ongoing with Quanterix

Increasing customer pull through

Consumable pull through from top 10
customers increased **3.5x since 2019**

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

Simoa biomarkers continue to enable therapy approvals



Failed Phase 3 study

108-subject study:
Clinical measures alone did not show benefit

OCT 2021

Quanterix NfL biomarker

provides evidence to show reduction in plasma neurofilament light

NFL

FDA accelerated approval

NfL provides enough basis for a conditional approval

APR 2023

Establishing Quanterix biomarker as the non-invasive proxy for brain health driving the neuro decade

Growing trials and approved therapies

2 drugs* approved

Academic Focus
~**2500** publications

Last 20 years

4 drugs* approved

All **4** fueled by **Quanterix**

- ✓ Revenue crossing **\$100M**
- ✓ **>45** active projects

Quanterix
Last 3 years

>120 biopharma CNS clinical-stage programs of which **24** programs focused specifically on AD


Long runway of continued double-digit growth in research

2024 and beyond



Multiple ongoing studies in Neurology

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



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

Alzheimer's - 1 in 9 people aged 65 and older in the US has it



>55M globally
living with Alzheimer's

By 2050 projected to
rise to 139M



Every 3 seconds
someone in the world
develops Alzheimer's



US\$1.3 trillion
global societal cost of
dementia in 2019



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



Detecting early with Simoa will impact quality of life

**Meets NIA-AA
Revised Criteria ...**

p-Tau 217

for Diagnosis and Staging
of Alzheimer's Disease

**... to address a
significant need**

75% people

not diagnosed
globally

Offering a test for all patients - LucentAD p-Tau 217

Alzheimer's Disease Continuum

p-Tau 217

EARLY STAGE

LATE STAGE

 Lucent Diagnostics

- ✓ 93% Accuracy
- ✓ 98% Specificity
- ✓ 90% Sensitivity



Earliest Stage




Scalable workflow




Highly Accurate

Only test meeting the NIA-AA criteria for diagnosing and staging of AD

Quanterix best positioned for Alzheimer's detection today

 **Quanterix**
Enabled DX lab

 **Non-Quanterix**
DX lab

**Quantitative range
in clinical cohorts**

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

Cost of test

Scalable platform

**Multiplex capabilities
(for differentiated DX)**

Simoa®

High

Yes

Low

<\$400/test

Yes

Yes

Mass Spectrometry

High

Yes

High

>\$1200/test

No

Limited

Chemiluminescence

Limited

No

Low

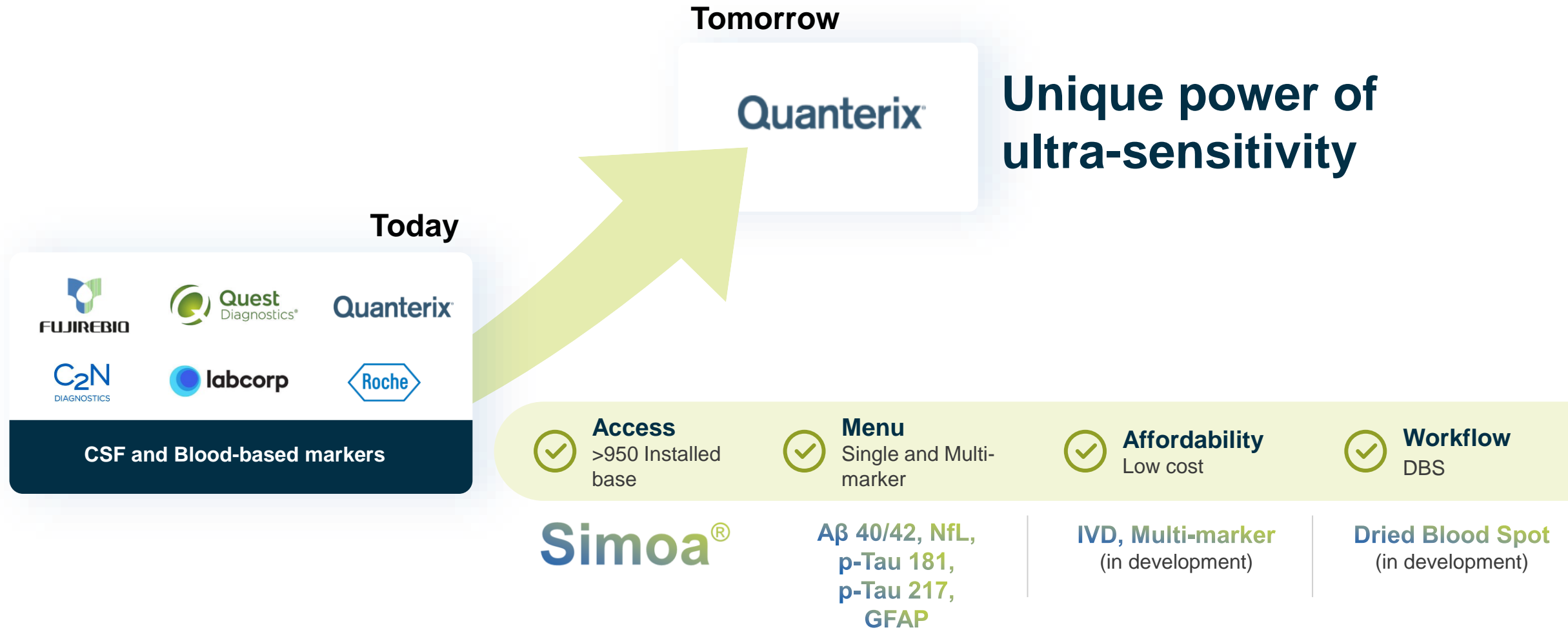
<\$400/test

Yes

No

**Simoa technology
uniquely suited for
building global
Diagnostics testing
infrastructure**

And uniquely set to lead the decade of Neuro



With a clear focus to maximize access to patients

Allocating capital

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

to execute commercial adoption

Reimbursement:
Submitted for CPT code (expected in 2024)

FDA:
Breakthrough designation received for 2
(expected 3rd in 2024)

Studies:
Four studies underway with additional
planned in 2024

Quanterix[®]

A diversified portfolio that is geared to **beat industry headwinds** with recent **transformation** and **multiple growth drivers**

Diversified portfolio

Mitigates macro headwinds in Life Science tools space by providing testing services as alternative

Indexed to well funded pharma

Several mid to late-stage programs continue to be funded providing recurring growth

Growing trials

While Dx ramps, continued growth in neuro trials (>124 CNS programs¹) will bolster research business for years to come

Built to lead neuro decade

Strong Results

Double digit top-line growth¹

Non-GAAP GM² of appr. 50%

Improvement of 1200 bps

Strong Balance Sheet

>\$300M¹ in net cash

Annual cash burn¹ down by >3x

Operational Leverage

Lab Services capability

>0.75M tests per year
(improvement of 50%)

Manufacturing capacity

>4M tests per year
(improvement of 300%)

Innovation rate

>20 assays per year
(development times reduced from >18 months to <6 months)

Clear Strategy

Research

Power the neuro decade with unique biomarkers

Diagnostics

Establish and grow the AD diagnostics business

1. Preliminary unaudited full year results for 2023.

2. The Company is guiding full year 2023 GAAP gross margin percentage to be in the high 50's, and non-GAAP gross margin percentage to be approximately 50%. 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.