

Quanterix

Powering a Revolution in Healthcare



```
struct group_info Init_groups = { .usage = ATOMIC_INIT(2) };  
struct group_info Init_groups = { .usage = ATOMIC_INIT(2) };  
struct group_info *groups_alloc(int gidsetsize){  
    struct group_info *groups_alloc(int gidsetsize){
```

Q1 2018 Quanterix Earnings Conference Call

- Q1 2018 Highlights: Kevin Hrusovsky Chairman & CEO - Slides
- Financial Results: Joe Driscoll CFO – Script
- Update and Q&A: Kevin Hrusovsky

Safe Harbor Statement



This presentation and the accompanying oral commentary contain “forward-looking” statements within the meaning of the Private Securities Litigation Reform Act of 1995, that are based on our beliefs and assumptions and on information available to us as of the date of this presentation. Forward-looking statements include statements that are not historical facts and can be identified by terms such as “anticipate,” “believe,” “continue,” “could,” “estimate,” “expect,” “intend,” “may,” “might,” “objective,” “ongoing,” “plan,” “predict,” “project,” “potential,” “should,” “will,” or “would,” or the negative of these terms, or other comparable terminology intended to identify statements about the future. We may not actually achieve the plans, intentions or expectations disclosed in our forward-looking statements. Forward-looking statements involve known and unknown risks, uncertainties, assumptions and other factors that may cause our actual results, performance or achievements to be materially different from any future results, performance or achievements expressed or implied by the forward-looking statements. The risks and uncertainties that we face are described in our most recent filings with the Securities and Exchange Commission. Except as required by law, we assume no obligation to update these forward-looking statements publicly, even if new information becomes available in the future.



Business

- Strong demand for SR-X platform, 1st. commercial quarter
- Record level of quarterly consumable revenue & instrument placements
- Closed Aushon Biosystems acquisition
- >100 Neuro publications, >215 in total, 500+ phase I-III trials, 18 of 20 top biopharmas
- Favorable macro trends for Biomarkers
- Finalist for FNIH Biomarker Consortium
- Strong advances in MS biomarkers
- Sales Team expansion for Pharma Services

Financial

- Revenue \$7.5M, +41% vs. PY Q1
- High margin business segments grew 60%, led by Neurology & now represent nearly 60% of total revenue
- Installed base continues to grow productively with avg annual consumable / inst exceeding \$50k/yr target (after six months).
- Pharma & Academics, +48% & 33% respectively
- Cost basis impacted by Aushon
- Expansion ongoing: HC 157, +31 vs. YE



Timing	Value Drivers and Opportunity
2018	CLIA Lab, Facilities, Menu-200 assays, Leadership / Team
2019	+ Pharma Services - Accelerator
2020	+ New Instruments and CDx - LDT National Coverage Decision
Longer-term	+ IVD and POC Entry

SR-X Interest Continues to Expand



Rapid Growth with Compelling Consumable Pull-through

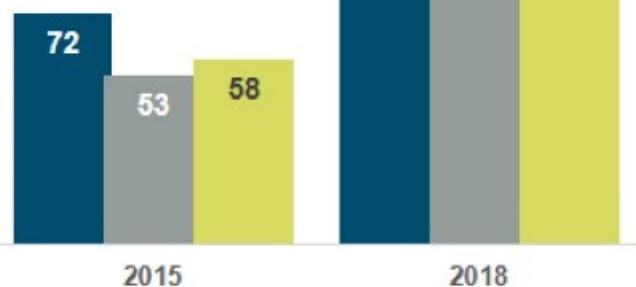


Portfolio Growth

(\$ in millions)

- Instruments
- Publications
- Markers

(cumulative)



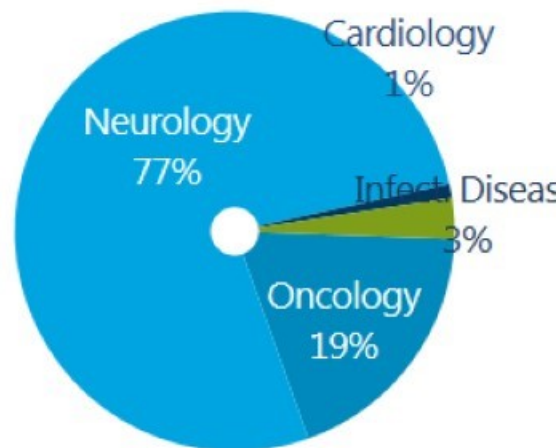
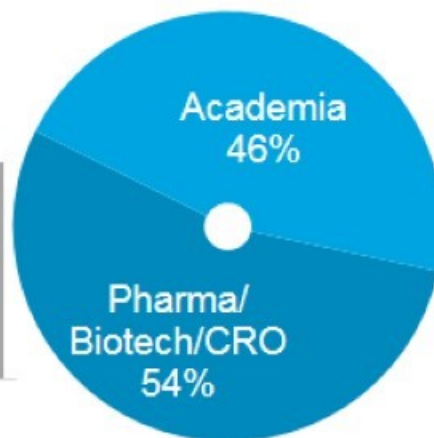
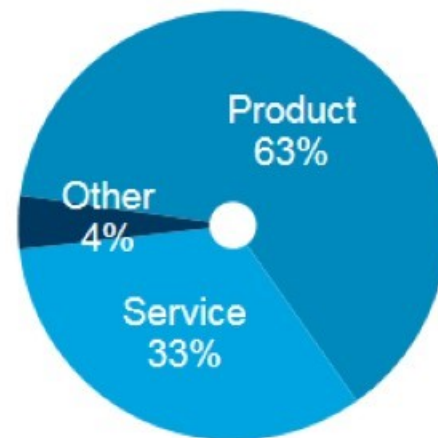
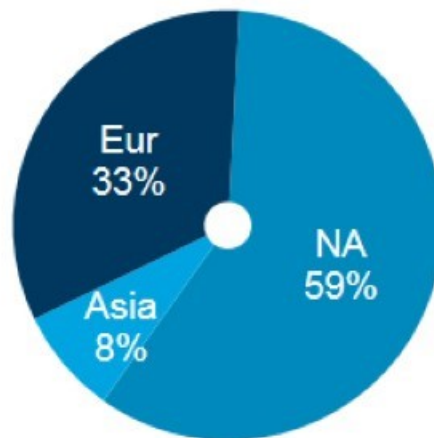
Grew from 0 instruments Jan'14

Revenue

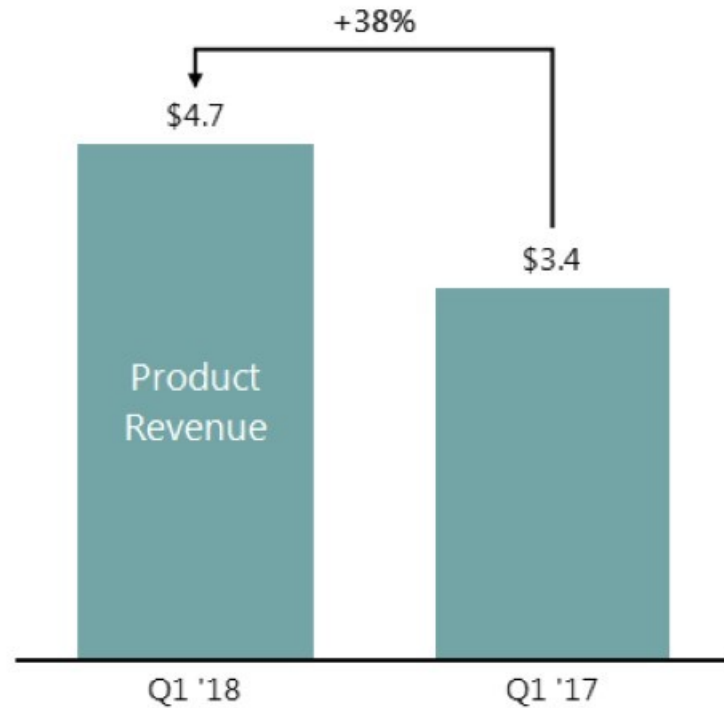
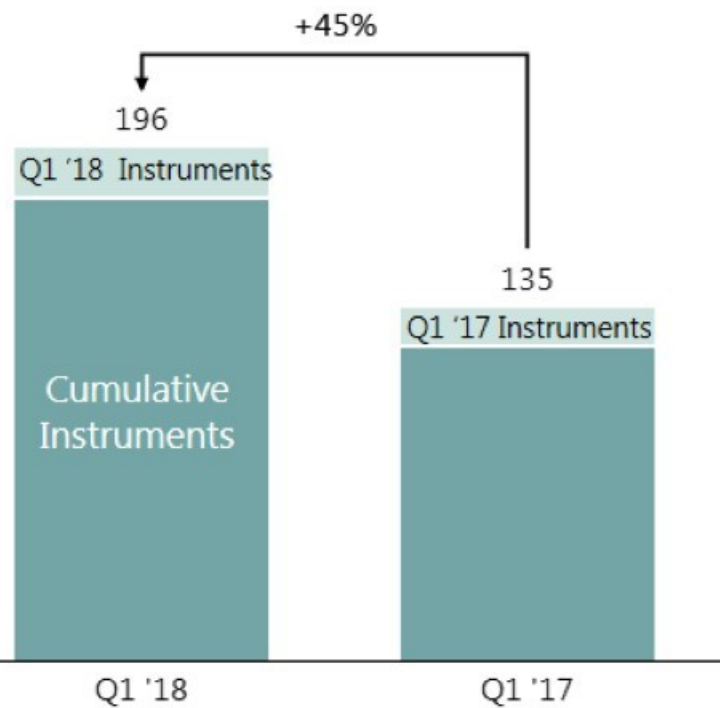


■ = Q1 Revenue

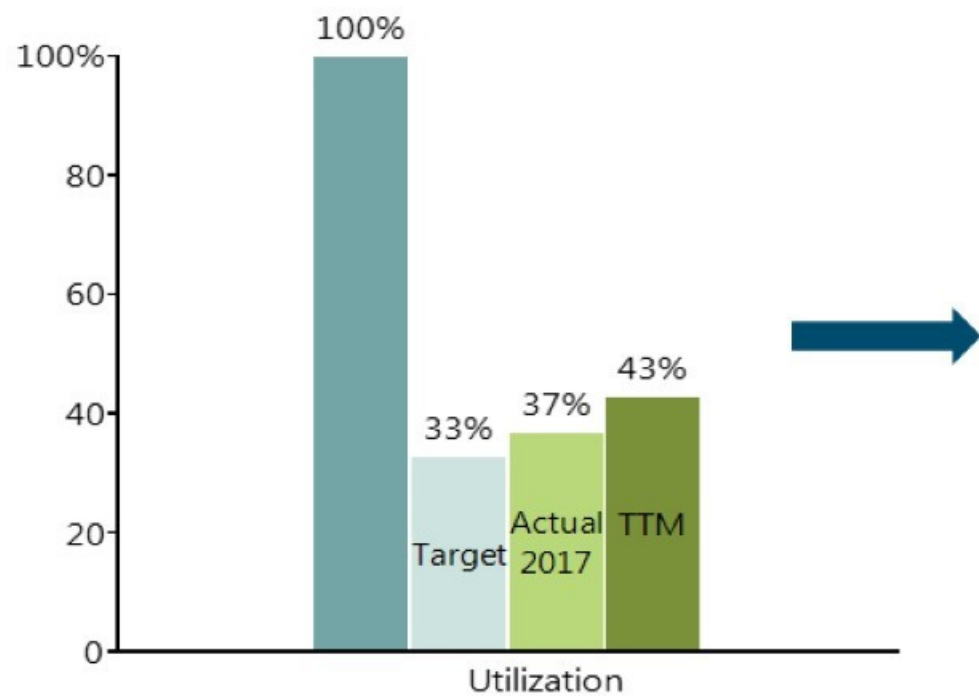
Q1 '18 Revenue



Strong Growth Across All Segments

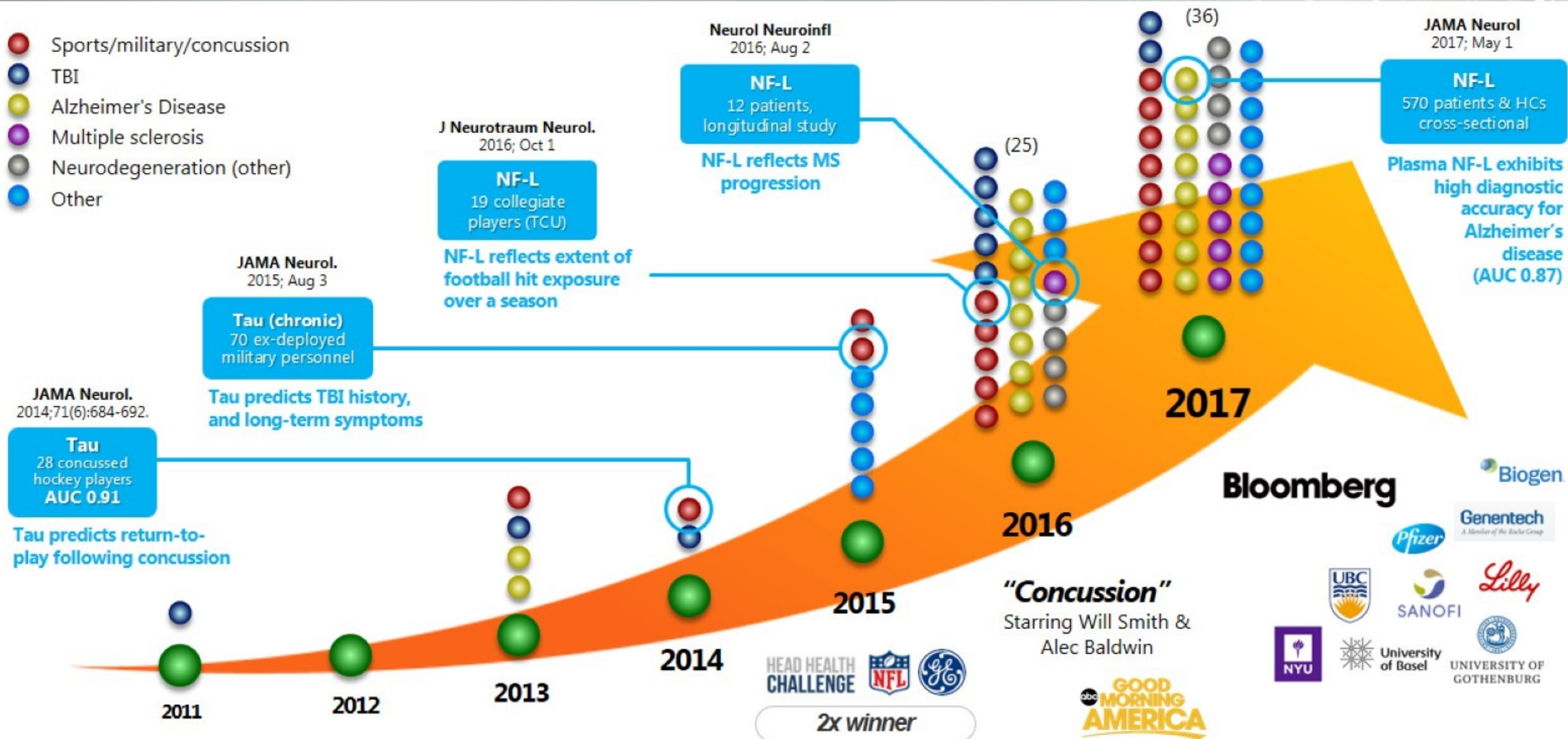


High Margin Consumable Revenue Continues to Scale



Consumable Growth:
Q1 '18 vs. Q1 '17 65%
FY '17 vs. FY '16 70%

Neurology Publications – Catalyze Disruption



Business-Favorable Changes in IVD Regulatory Climate



Early Alzheimer's Disease: Developing Drugs for Treatment Guidance for Industry



FDA initiative...

FDA recognizes the criticality of ID'ing *pre-clinical* patients for drug trials using biomarkers.

What it means for QTRX...

Encourages biopharma to identify/incorporate promising biomarkers in trial design

Contains Nonbinding Recommendations
Draft - Not for Implementation

Breakthrough Devices Program Draft Guidance for Industry and Food and Drug Administration Staff



FDA to expedite development, assessment, and review of medical devices addressing treatment or diagnosis of life-threatening or irreversibly debilitating conditions.

Expedited review of IVD tests; more probable Class II path. Example: Banyan TBI test (6 mo to clearance).

FEDERAL REGISTER The Daily Journal of the United States Government

Program for Parallel Review of Medical Devices

A Notice by the Centers for Medicare & Medicaid Services and the Food and Drug Administration on 10/24/2016

Leroy A. Richardson,
Chief, Information Collection Review Office,
Office of Scientific Integrity, Office of the
Associate Director for Science, Office of the
Director, Centers for Disease Control and
Prevention.
FR Doc. 2016-21001 Filed 10-25-16; 8:41 am
BILLING CODE 4150-10-P

DEPARTMENT OF HEALTH AND
HUMAN SERVICES

Centers for Medicare & Medicaid
Services

[CMS-3180-H4]

Food and Drug Administration

[Docket No. FDA-2010-N-0308]

Program for Parallel Review of Medical
Devices

AGENCY: Food and Drug Administration;
Centers for Medicare & Medicaid
Services, HHS.

ACTION: Notice.

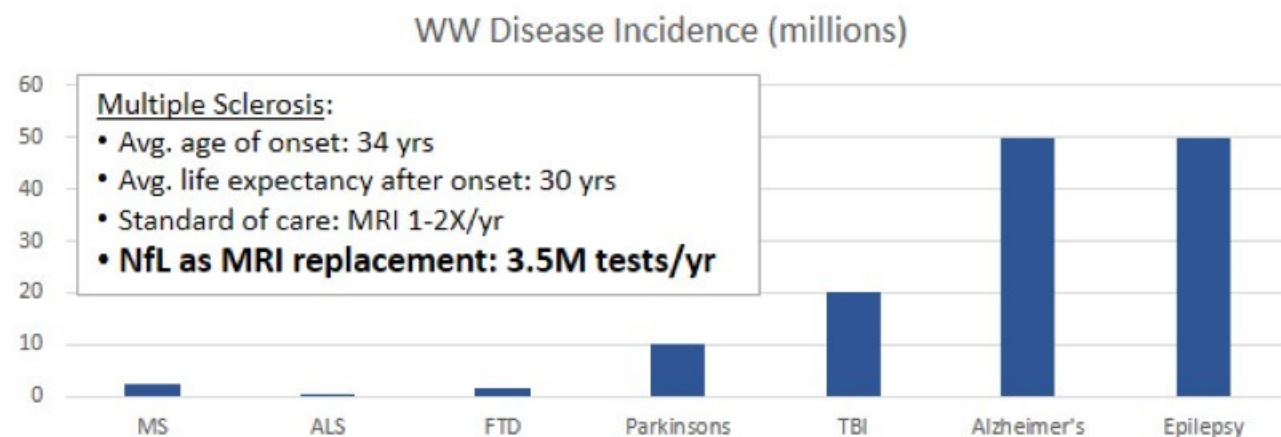
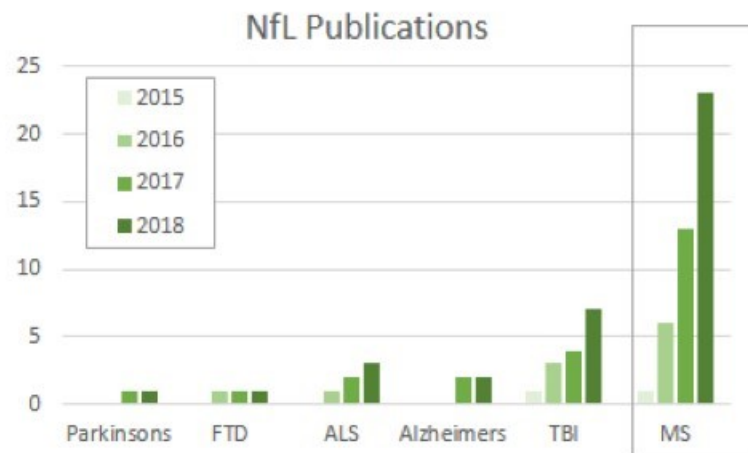
SUMMARY: The Food and Drug Administration (FDA) and the Centers for Medicare & Medicaid Services (CMS) (the Agencies) are informing the public that the Parallel Review of medical devices pilot program will be fully implemented and extended indefinitely. The Agencies are soliciting nominations from manufacturers of innovative medical devices to participate in the "Program for Parallel Review of Medical Devices." The Parallel Review program is a collaborative effort that is intended to reduce the time between FDA marketing approval or FDA's granting of a *de novo* request and Medicare coverage decisions through CMS's National Coverage Determination (NCD).



FDA teaming up with CMS to provide parallel review of medical device approval and Medicare coverage.

Potential to secure reimbursement determination simultaneous with FDA clearance

Emerging Clinical Biomarker: Neurofilament Light (NfL)



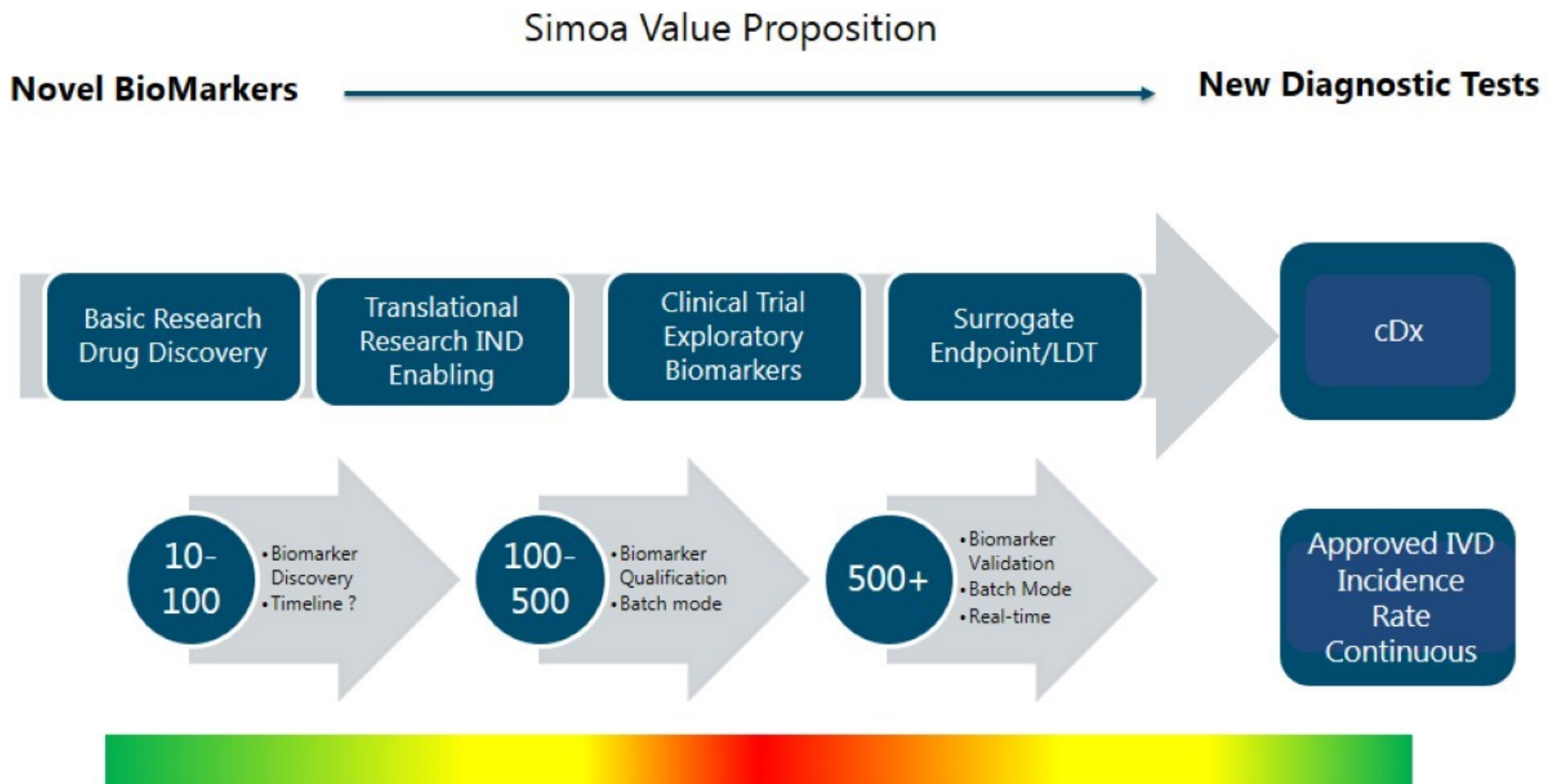
Studies confirm NfL clinical utility:

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

Majority of published data obtained with Simoa NfL

Clinical Validation of NfL for MS is a Key Beachhead

Quanterix Biomarker Support for Clinical Trials



Representative Accelerator Projects



Sponsor	Indication	Phase I	Phase II	Phase III	Current \$	CDx Potential
Pharma 1	Mood Disorders				\$2.2M	✓
Pharma 2	Asthma				\$900M	✓
Biotech 1	MS				\$500M	
Pharma 1	Alzheimer's				\$450M	
Pharma 3	MS				\$320M	✓
Pharma 1	Alzheimer's				\$250M	
Pharma 4	MS				\$250M	
Biotech 1	MS				\$200M	✓
Biotech 2	Asthma			PD	\$200M	
Biotech 3	Crohn's				\$100M	✓



Joe Driscoll CFO provides Q1 2018 Financial Results - Script

PREVENT · DETECT · TREAT

Save the Date
3rd Annual
Powering Precision Health Summit 2018
December 11-12, 2018
Boston Marriot Newton
Targeting 1000 + attendees, 65 Speakers, 25 Sponsors

POWERING PRECISION HEALTH

“If we start the PPH movement now, by 2035, healthcare will be 40% lower in cost, 60% more accessible to average citizens, and we will live 8 years longer productive lives.”

Kevin Hrusovsky

Founder & Chair, Powering Precision Health Summit

606

attendees

54

luminary speakers

18

corporate sponsors

Revolutionary opportunities to detect earlier and improve treatment of:

- Pancreatic Cancer
- Breast Cancer
- Prostate Cancer
- Lung Cancer
- Alzheimer's
- CTE
- Concussions
- Diabetes
- Multiple Sclerosis
- Parkinson
- HIV
- C-Difficile

Speakers from:

Harvard Medical, MD Anderson Cancer Center, Merck, Eli Lilly, NYU, UPenn, NIH, DanaFarber, Yale Medical, Biogen, Cedars Sinai, Myriad RBM, Boston Univ.

Massachusetts Governor Issues Proclamation for Precision Health Awareness Week

12

PPH news articles

402

tweets about PPH

154

live streams on Periscope

133,000,000

social media impressions

Quanterix

Powering a Revolution in Healthcare



```
struct group_info Init_groups = { .usage = ATOMIC_INIT(2) };  
struct group_info Init_groups = { .usage = ATOMIC_INIT(2) };  
struct group_info *groups_alloc(int gidsetsize){  
    struct group_info *groups_alloc(int gidsetsize){
```

2018 Update and Q&A: Kevin Hrusovsky



Poised to Disrupt Healthcare and Create Significant Value



¹ \$30bn subject to receipt of regulatory approvals or clearances, which the company has not applied for to date.