

Product Number:	103345
Lot Number:	501826
Expiration:	8-May-2020
Platform(s):	HD-1 & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103338	916805
Detector Reagent	103339	916804
SBG Reagent	103340	912802
Sample Diluent	103341	912803
Calibrators	103342	910101
Controls	103343-103344	910102
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex B SR-X Kit Instructions (KI-0123).

Release Materials

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	1.89	1.28-2.36
Control 2	910102	94.8	73.8-111
Panel 1	MF-101617-01	1.57	1.06-1.75
Panel 2	MF-101617-02	37.6	27.4-41.1
Panel 3	MF-101617-03	151	117-175

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	5.38	4.59-6.88
Control 2	910102	545	453-679
Panel 1	MF-101617-01	4.00	2.81-4.60
Panel 2	MF-101617-02	38.0	22.3-40.8
Panel 3	MF-101617-03	373	313-469

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	104	80.4-121
Control 2	910102	11343	8772-13157
Panel 1	MF-101617-01	49.2	26.7-53.1
Panel 2	MF-101617-02	165	128-192
Panel 3	MF-101617-03	2452	2118-3177

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	109	85.2-135
Control 2	910102	10458	8319-12478
Panel 1	MF-101617-01	64.4	42.2-84.7
Panel 2	MF-101617-02	462	235-636
Panel 3	MF-101617-03	5152	4273-6409

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Patrizia Stadler Assoc. Director, QC

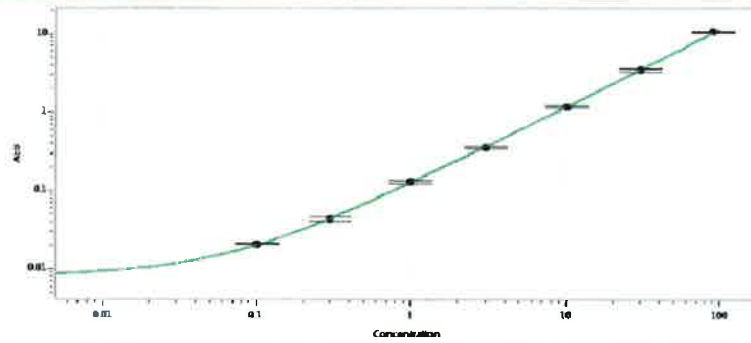
P. Stadler 20 Jun 2019

Name

Title

Signature/Date

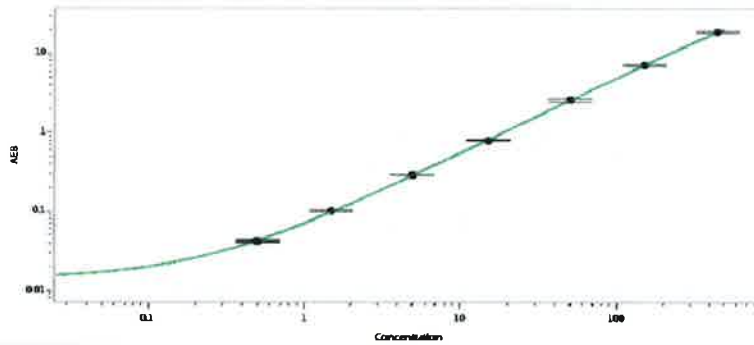
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.086
C	0.291
D	0.974
E	2.76
F	9.24
G	27.0
H	85.5

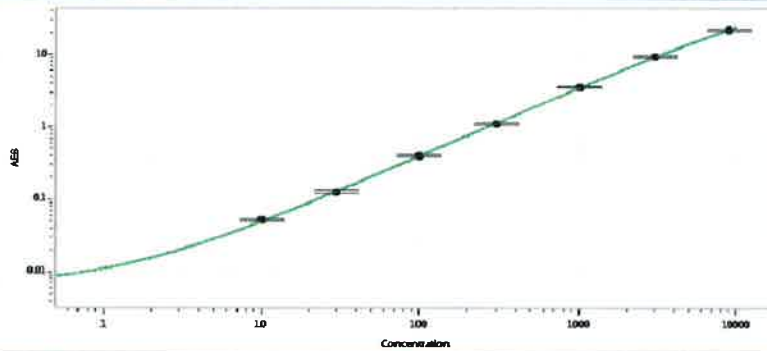
NF-light® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.466
C	1.49
D	5.02
E	14.9
F	51.9
G	149
H	451

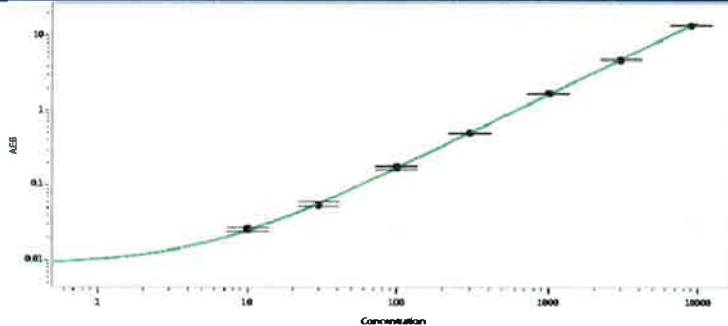
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	9.30
C	26.6
D	94.0
E	271
F	984
G	2800
H	8563

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	9.93
C	27.7
D	94.5
E	276
F	955
G	2753
H	8387