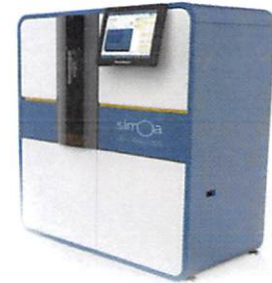


Product Number:	103345
Lot Number:	501712
Expiration:	12-Nov-2019
Platform(s):	HD-1 and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103338	902516
Detector Reagent	103339	902512
SBG Reagent	103340	831602
Sample Diluent	103341	831603
Calibrators	103342	910101
Controls	103343-103344	910102
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	1.81	1.24-2.37
Control 2	910102	91.4	73.4-110
Panel 1	MF-101617-01	1.36	1.06-1.69
Panel 2	MF-101617-02	31.0	26.8-40.2
Panel 3	MF-101617-03	151	116-174

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	5.84	4.63-6.95
Control 2	910102	550	456-683
Panel 1	MF-101617-01	3.95	2.75-4.56
Panel 2	MF-101617-02	25.6	23.9-35.9
Panel 3	MF-101617-03	347	315-472

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	103	79.3-121
Control 2	910102	10409	8721-13082
Panel 1	MF-101617-01	42.1	29.8-46.5
Panel 2	MF-101617-02	153	127-191
Panel 3	MF-101617-03	2587	2145-3218

UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910102	114	84.0-136
Control 2	910102	9994	8311-12466
Panel 1	MF-101617-01	67.3	41.0-85.5
Panel 2	MF-101617-02	342	211-668
Panel 3	MF-101617-03	4839	4299-6448

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

Review/Approval

Patrizia Stadler QC Assoc. Director

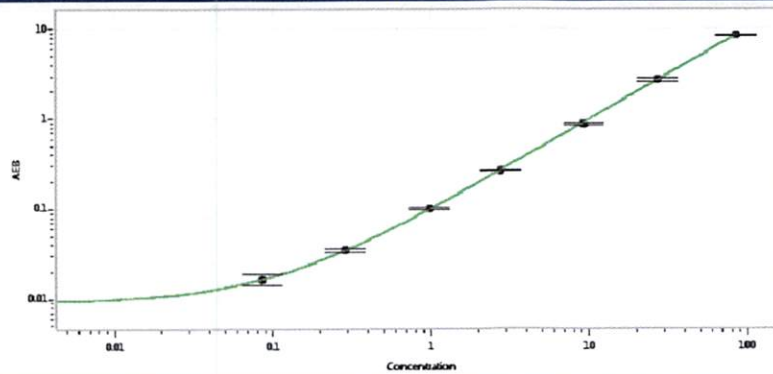
Name

Title

Signature/Date

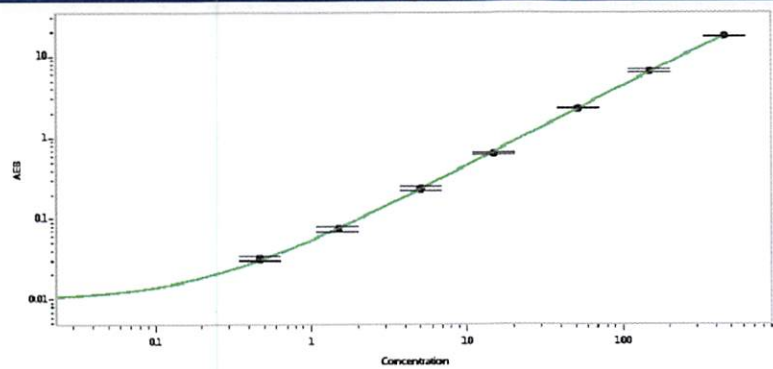
P. Stadler 29 Apr 2019

Tau Calibration Curve



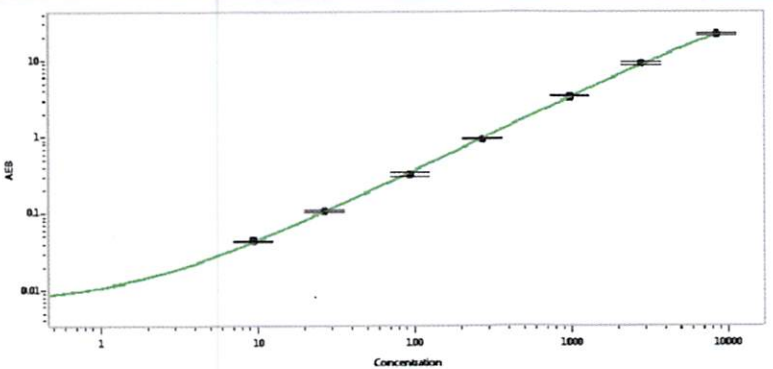
Calibrator Levels (pg/mL)	
A	0.000
B	0.086
C	0.291
D	0.974
E	2.76
F	9.24
G	27.0
H	85.5

NF-light® Calibration Curve



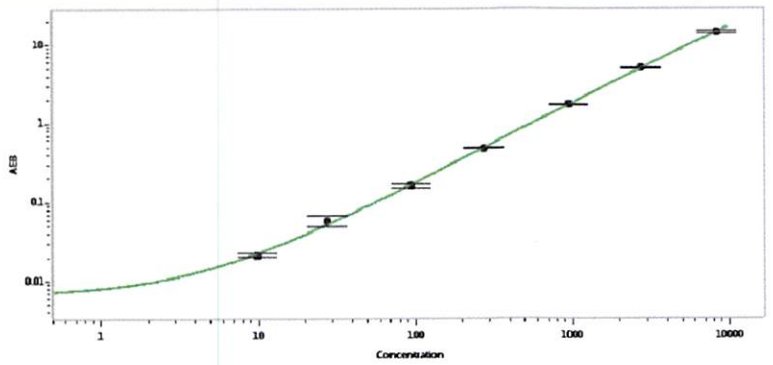
Calibrator Levels (pg/mL)	
A	0.000
B	0.466
C	1.49
D	5.02
E	14.9
F	51.9
G	149
H	451

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