

Product Number:	102153
Lot Number:	501687
Expiration:	19-Feb-2020
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	102141	908703
Detector Reagent	102143	908704
SBG Reagent	102145	908604
Sample Diluent	102147	905009
Calibrators	102150	905901
Controls	102151-102152	905902
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	905902	1.85	1.57-2.36
Control 2	905902	90.0	76.3-114
Panel 1	MF-101617-01	1.09	0.710-1.65
Panel 2	MF-101617-02	32.9	22.7-44.9
Panel 3	MF-101617-03	152	93.3-204

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	905902	5.32	3.95-5.92
Control 2	905902	479	392-589
Panel 1	MF-101617-01	4.23	3.08-6.35
Panel 2	MF-101617-02	38.4	16.8-50.2
Panel 3	MF-101617-03	386	232-548

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	905902	23.0	17.9-26.8
Control 2	905902	1116	901-1352
Panel 1	MF-101617-01	48.2	40.6-62.2
Panel 2	MF-101617-02	195	110-229
Panel 3	MF-101617-03	2640	2138-3313

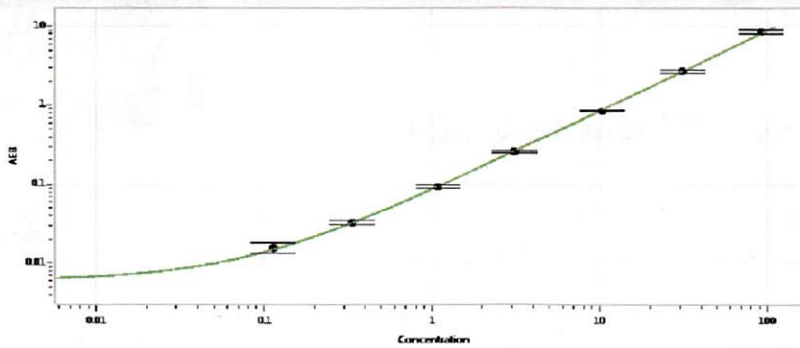
UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	905902	94.6	69.4-126
Control 2	905902	8854	7226-10838
Panel 1	MF-101617-01	63.7	39.0-89.1
Panel 2	MF-101617-02	445	308-473
Panel 3	MF-101617-03	5044	4253-6380

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

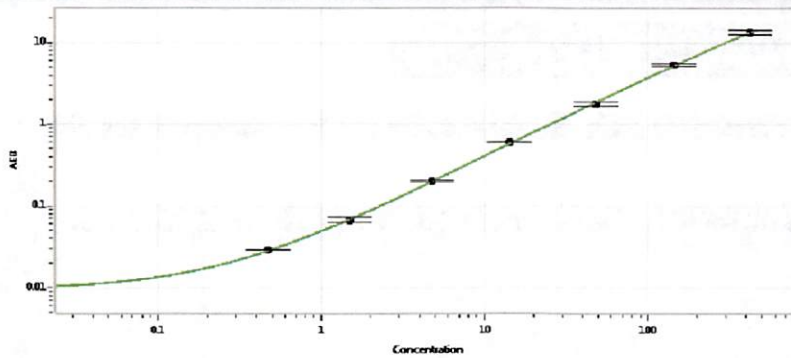
Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	09 April 2019
	Name	Title	Signature/Date	

Tau Calibration Curve



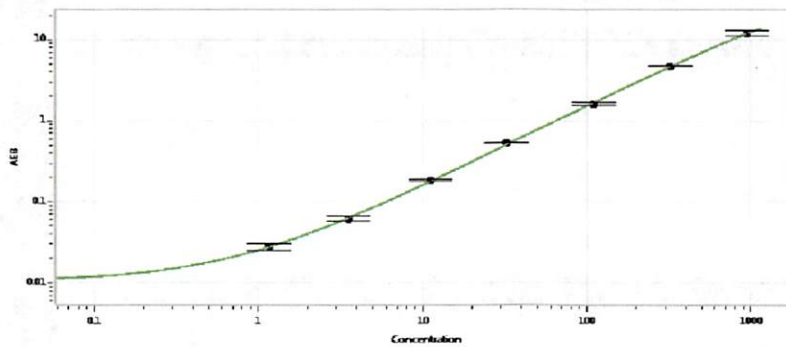
Calibrator Levels (pg/mL)	
A	0.000
B	0.113
C	0.332
D	1.08
E	3.12
F	10.3
G	31.3
H	91.6

NF-light® Calibration Curve



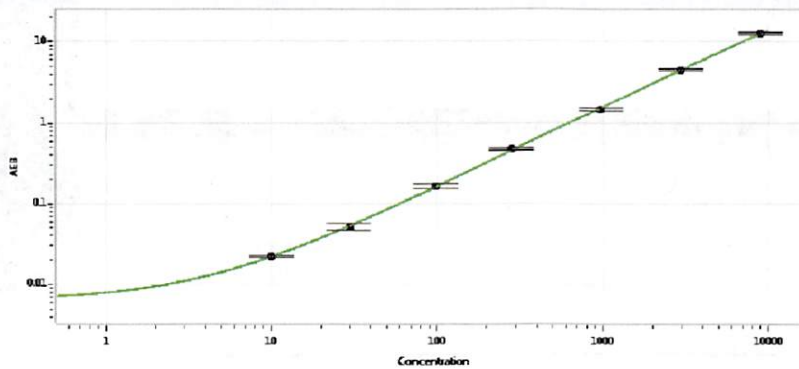
Calibrator Levels (pg/mL)	
A	0.000
B	0.470
C	1.50
D	4.80
E	14.3
F	48.0
G	144
H	426

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.16
C	3.54
D	11.1
E	32.6
F	111
G	323
H	946

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	10.0
C	29.6
D	97.9
E	285
F	980
G	2940
H	8998