

Product Number:	102153
Lot Number:	501762
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	909505
SBG Reagent	102145	908604
Sample Diluent	102147	905009
Calibrators	102150	905901
Controls	102151-102152	905902
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	2.10	1.57-2.36
Control 2	905902	94.8	76.3-114
Panel 1	MF-101617-01	1.15	0.710-1.65
Panel 2	MF-101617-02	34.1	22.7-44.9
Panel 3	MF-101617-03	150	93.3-204

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	905902	5.45	3.95-5.92
Control 2	905902	513	392-589
Panel 1	MF-101617-01	4.65	3.08-6.35
Panel 2	MF-101617-02	39.0	16.8-50.2
Panel 3	MF-101617-03	398	232-548

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	905902	25.5	17.9-26.8
Control 2	905902	1147	901-1352
Panel 1	MF-101617-01	51.4	40.6-62.2
Panel 2	MF-101617-02	197	110-229
Panel 3	MF-101617-03	2617	2138-3313

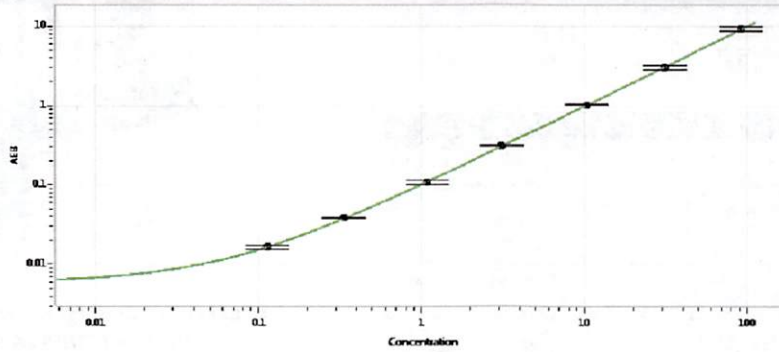
**UCHL1®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	905902	92.6	69.4-126
Control 2	905902	9199	7226-10838
Panel 1	MF-101617-01	61.6	39.0-89.1
Panel 2	MF-101617-02	423	308-473
Panel 3	MF-101617-03	5228	4253-6380

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

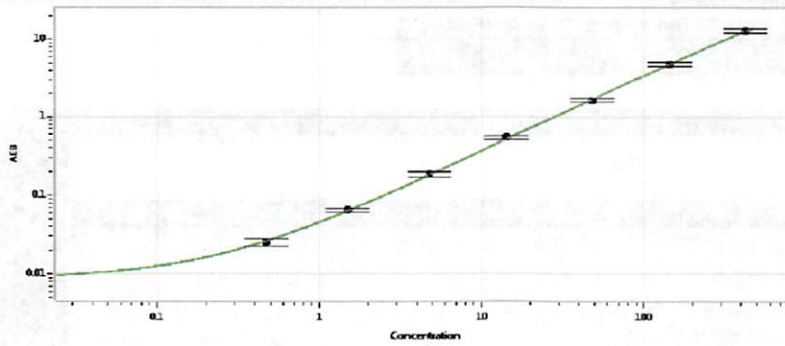
Review/Approval	Patrizia Stadler QC Assoc. Director	<i>P. Stodler</i>	10 Apr. 1 2019
	Name Title	Signature/Date	

**Tau Calibration Curve**



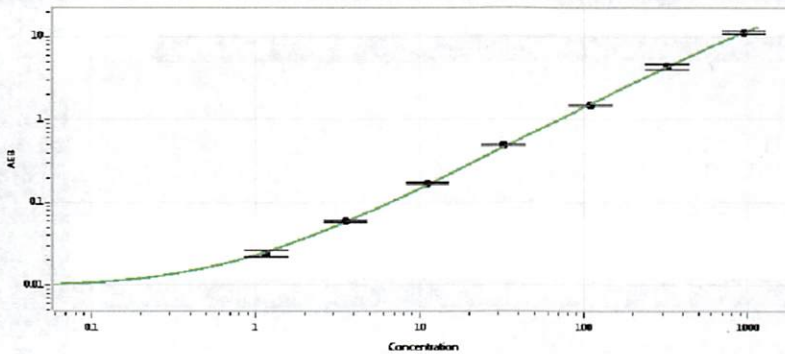
Calibrator Levels (pg/mL)	
A	0.000
B	0.113
C	0.332
D	1.08
E	3.12
F	10.3
G	31.3
H	91.6

**NF-light® Calibration Curve**



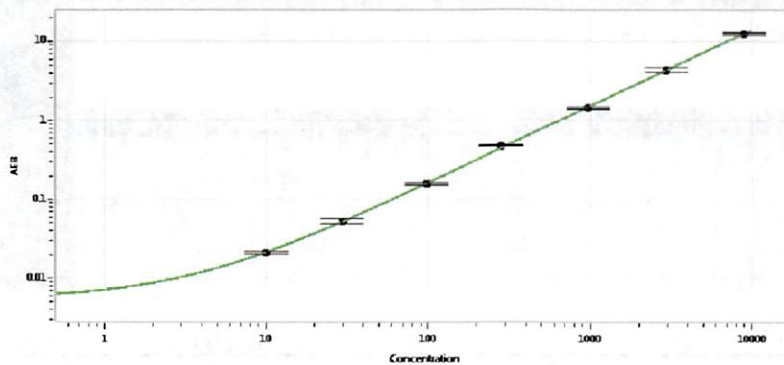
Calibrator Levels (pg/mL)	
A	0.000
B	0.470
C	1.50
D	4.80
E	14.3
F	48.0
G	144
H	426

**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