

Product Number:	102153
Lot Number:	501583
Expiration:	10-Jan-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	902515
Detector Reagent	102143	902511
SBG Reagent	102145	901004
Sample Diluent	102147	901105
Calibrators	102150	827401
Controls	102151-102152	827402
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	827402	2.15	1.36-2.88
Control 2	827402	99.8	54.8-136
Panel 1	MF-101617-01	1.15	0.710-1.65
Panel 2	MF-101617-02	34.8	22.7-44.9
Panel 3	MF-101617-03	134	93.3-204

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	5.00	3.78-7.36
Control 2	827402	514	267-719
Panel 1	MF-101617-01	4.23	3.08-6.35
Panel 2	MF-101617-02	29.8	16.8-50.2
Panel 3	MF-101617-03	348	232-548

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	22.0	17.0-25.5
Control 2	827402	1060	629-1390
Panel 1	MF-101617-01	48.9	40.6-62.2
Panel 2	MF-101617-02	197	110-229
Panel 3	MF-101617-03	2624	2138-3313

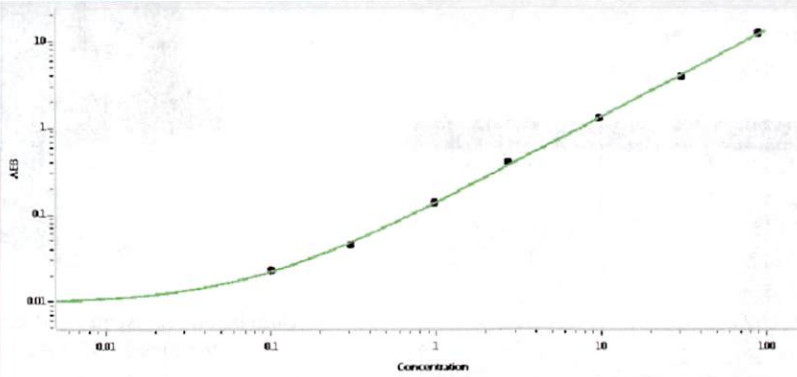
**UCHL1®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	121	89.2-151
Control 2	827402	11475	6829-14447
Panel 1	MF-101617-01	64.3	39.0-89.1
Panel 2	MF-101617-02	360	308-473
Panel 3	MF-101617-03	5255	4253-6380

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

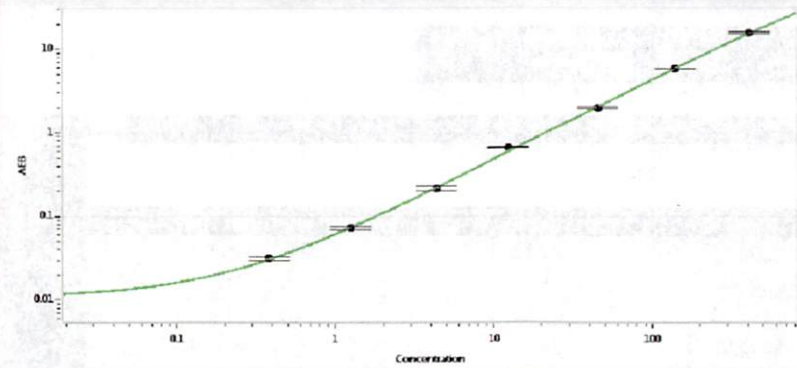
Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	30 Jan 2019
	Name	Title	Signature/Date	

**Tau Calibration Curve**



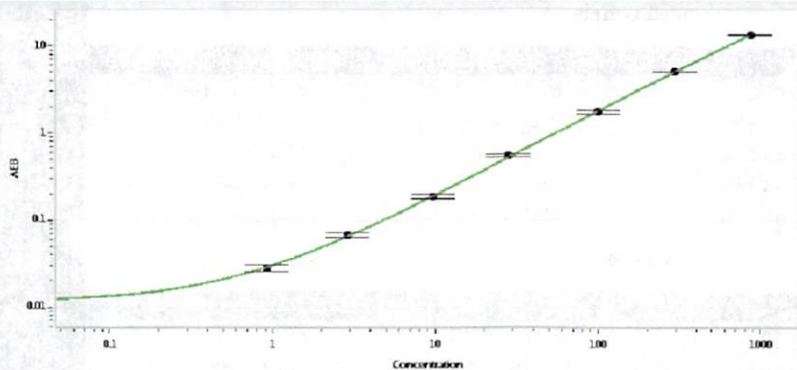
Calibrator Levels (pg/mL)	
A	0.000
B	0.100
C	0.302
D	0.986
E	2.73
F	9.80
G	30.3
H	89.8

**NF-light® Calibration Curve**



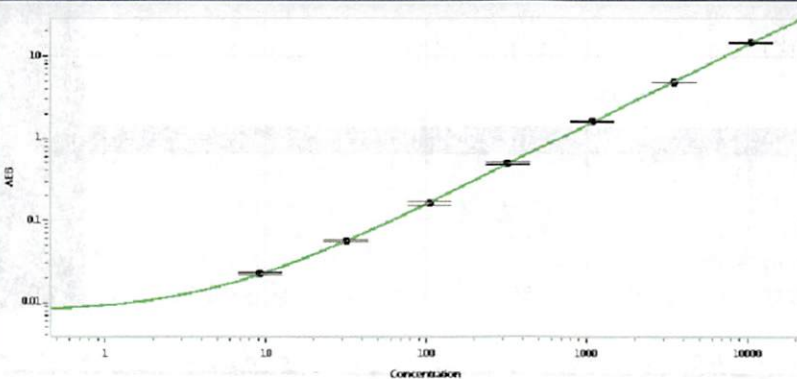
Calibrator Levels (pg/mL)	
A	0.000
B	0.380
C	1.26
D	4.36
E	12.4
F	45.3
G	139
H	408

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.923
C	2.91
D	9.73
E	28.2
F	99.8
G	298
H	866

**UCHL1® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	9.14
C	31.9
D	106
E	320
F	1095
G	3488
H	10606