

Product Number:	102153
Lot Number:	501380
Expiration:	10-Feb-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	102141	823404
Detector Reagent	102143	823309
SBG Reagent	102145	822201
Sample Diluent	102147	822202
Calibrators	102150	81710/821104
Controls	102151-102152	821105
SBG Diluent (SR-X)†	103158	N/A <sup>†</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	821105	2.08	1.71-2.65
Control 2	821105	109	85.3-128
Panel 1	MF-101617-01	1.17	0.710-1.65
Panel 2	MF-101617-02	32.6	22.7-44.9
Panel 3	MF-101617-03	146	93.3-204

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	821105	5.11	3.53-7.75
Control 2	821105	497	408-612
Panel 1	MF-101617-01	4.40	3.08-6.35
Panel 2	MF-101617-02	28.1	16.8-50.2
Panel 3	MF-101617-03	320	232-548

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	821105	21.1	17.5-26.5
Control 2	821105	1113	904-1356
Panel 1	MF-101617-01	60.4	40.6-62.2
Panel 2	MF-101617-02	167	110-229
Panel 3	MF-101617-03	2483	2138-3313

**UCHL1®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	821105	124	90.9-141
Control 2	821105	11037	9026-13540
Panel 1	MF-101617-01	58.3	39.0-89.1
Panel 2	MF-101617-02	371	308-473
Panel 3	MF-101617-03	5707	4253-6380

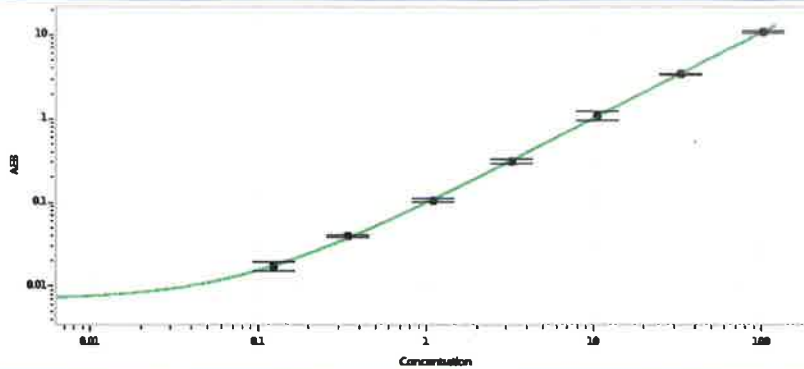
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Sivasankari Murugan QC Supervisor  
Name Title

*S. Murugan* 06 SEP 2018  
Signature/Date

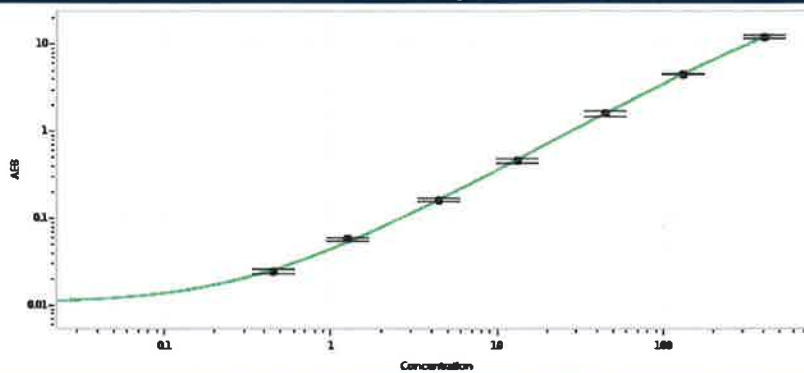
**Tau Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.124
C	0.341
D	1.10
E	3.24
F	10.5
G	33.0
H	103

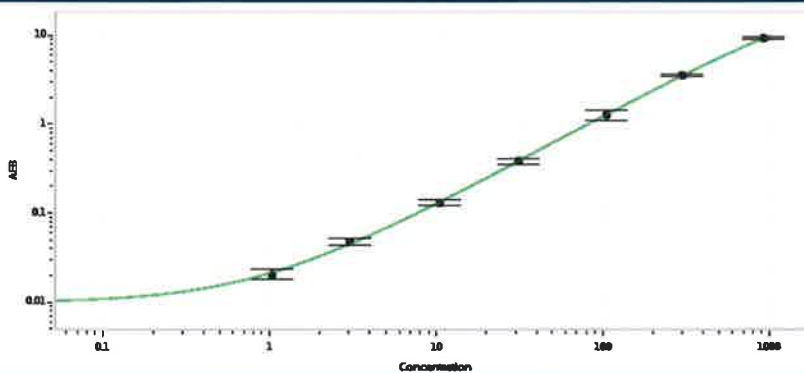
**NF-light® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.451
C	1.27
D	4.39
E	13.1
F	44.6
G	132
H	406

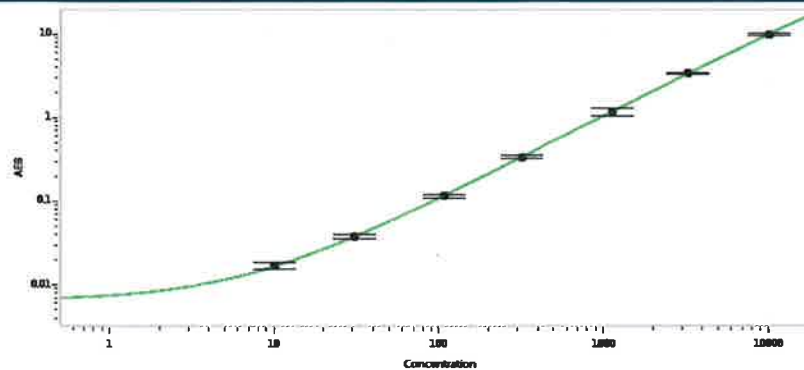
**GFAP® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	1.03
C	3.02
D	10.4
E	31.0
F	106
G	298
H	929

**UCHL1® Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	10.0
C	30.7
D	108
E	320
F	1133
G	3256
H	10168