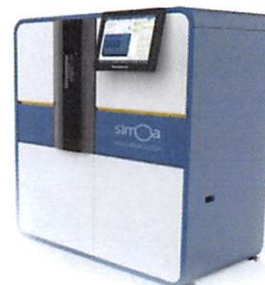


Product Number:	102153
Lot Number:	501443
Expiration:	10-Feb-2019
Platform(s):	HD-1 and SR-X



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	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	2.00	1.32-2.90
Control 2	827402	91.1	52.3-137
Panel 1	MF-101617-01	1.04	0.710-1.65
Panel 2	MF-101617-02	31.5	22.7-44.9
Panel 3	MF-101617-03	143	93.3-204

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	4.86	3.82-7.45
Control 2	827402	503	254-728
Panel 1	MF-101617-01	3.61	3.08-6.35
Panel 2	MF-101617-02	26.4	16.8-50.2
Panel 3	MF-101617-03	353	232-548

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	21.1	16.9-25.4
Control 2	827402	1056	606-1402
Panel 1	MF-101617-01	51.0	40.6-62.2
Panel 2	MF-101617-02	154	110-229
Panel 3	MF-101617-03	2617	2138-3313

**UCHL1®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	116	87.7-153
Control 2	827402	11045	6602-14489
Panel 1	MF-101617-01	59.6	39.0-89.1
Panel 2	MF-101617-02	362	308-473
Panel 3	MF-101617-03	5594	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

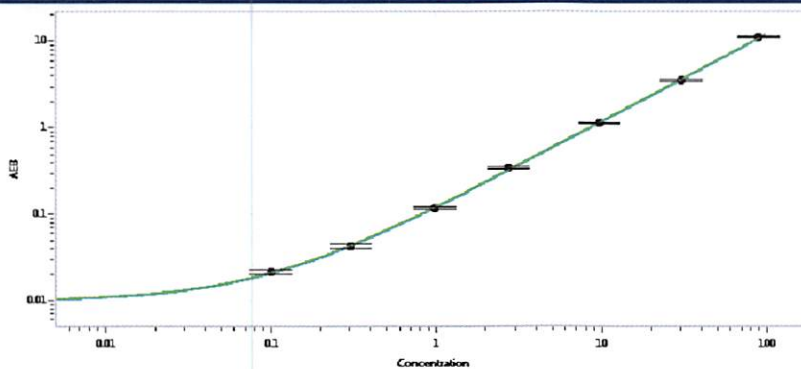
Name

Title

*[Signature]* 11 Oct 2018

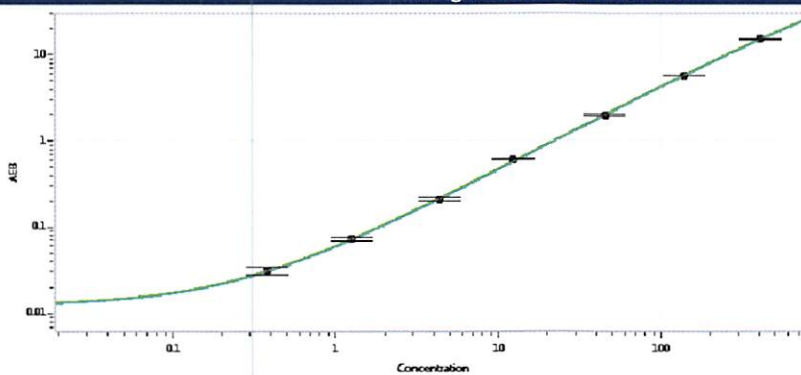
Signature/Date

**Tau Calibration Curve**



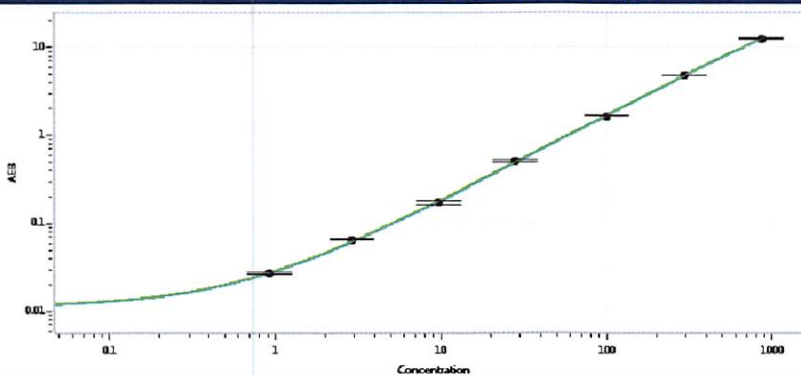
Calibrator Levels (pg/mL)	
A	0.000
B	0.100
C	0.302
D	0.986
E	2.73
F	9.80
G	30.3
H	89.8

**NF-light® Calibration Curve**



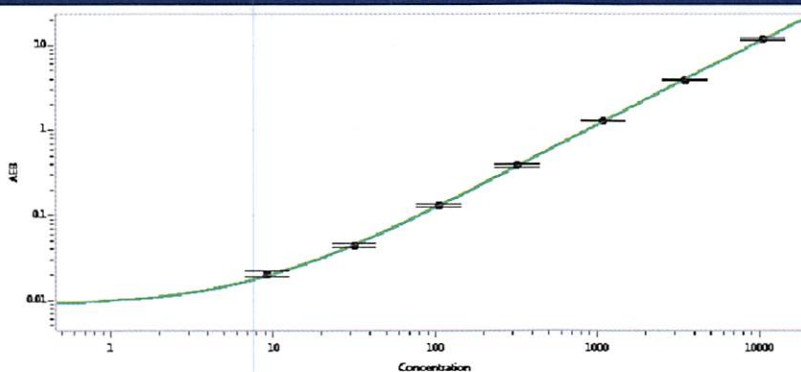
Calibrator Levels (pg/mL)	
A	0.000
B	0.380
C	1.26
D	4.36
E	12.4
F	45.3
G	139
H	408

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.923
C	2.91
D	9.73
E	28.2
F	100
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