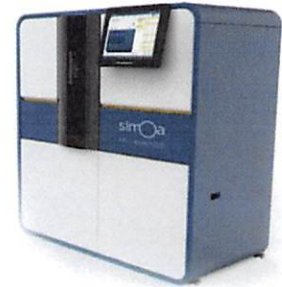


Product Number:	102153
Lot Number:	501428
Expiration:	21-Sep-2019
Platform(s):	HD-1 and SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102141	833908
Detector Reagent	102143	833907
SBG Reagent	102145	826411
Sample Diluent	102147	826502
Calibrators	102150	827401
Controls	102151-102152	827402
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	827402	2.03	1.32-2.90
Control 2	827402	96.6	52.3-137
Panel 1	MF-101617-01	1.07	0.710-1.65
Panel 2	MF-101617-02	30.8	22.7-44.9
Panel 3	MF-101617-03	142	93.3-204

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	5.34	3.82-7.45
Control 2	827402	496	254-728
Panel 1	MF-101617-01	4.09	3.08-6.35
Panel 2	MF-101617-02	30.8	16.8-50.2
Panel 3	MF-101617-03	370	232-548

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	22.5	16.9-25.4
Control 2	827402	995	606-1402
Panel 1	MF-101617-01	50.4	40.6-62.2
Panel 2	MF-101617-02	137	110-229
Panel 3	MF-101617-03	2240	2138-3313

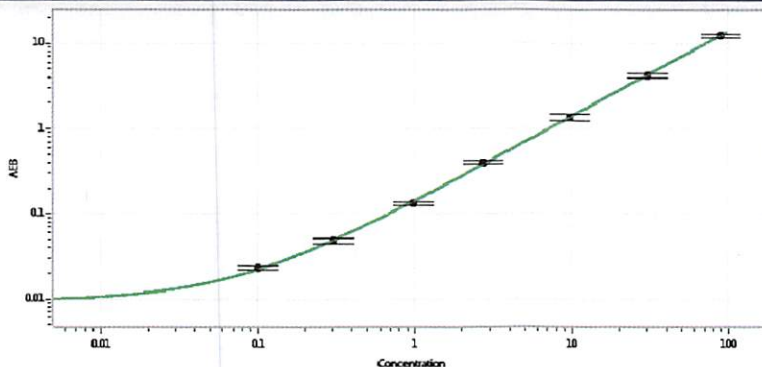
UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	827402	106	87.7-153
Control 2	827402	10899	6602-14489
Panel 1	MF-101617-01	48.7	39.0-89.1
Panel 2	MF-101617-02	298	308-473
Panel 3	MF-101617-03	4903	4253-6380

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

Review/Approval: Sivasankari Murugan (Name), QC Supervisor (Title), [Signature] 12DEC2018 (Signature/Date)

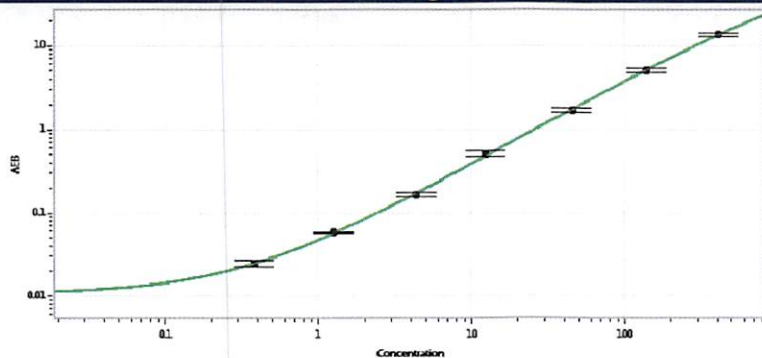
Tau Calibration Curve



**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

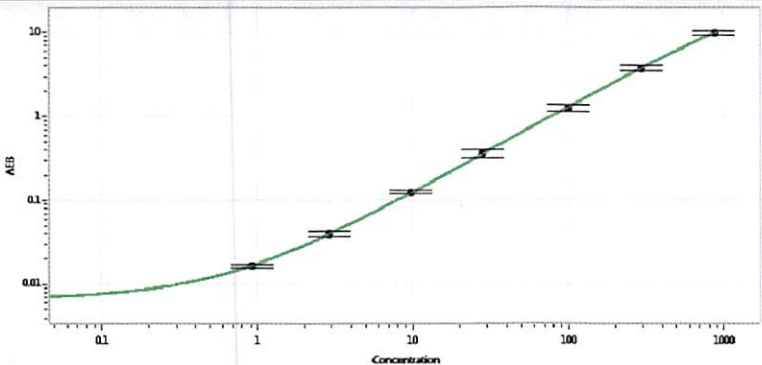
NF-light® Calibration Curve



**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

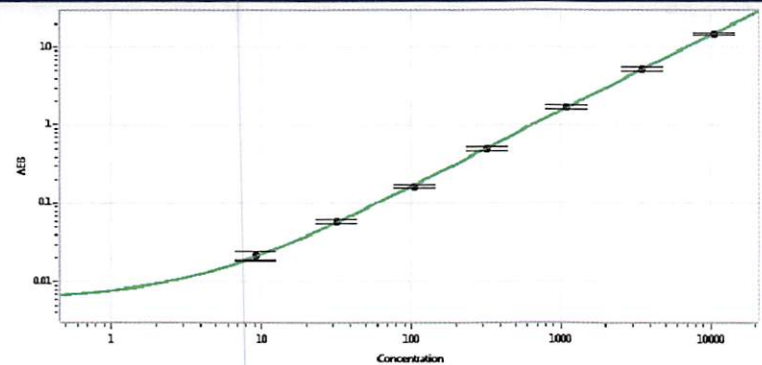
GFAP® Calibration Curve



**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