

Product Number:	102153
Lot Number:	501354
Expiration:	19-Dec-2018
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	820107
Detector Reagent	102143	820301
SBG Reagent	102145	81702
Sample Diluent	102147	81701
Calibrators	102150	81710/821104
Controls	102151-102152	821105
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	821105	2.32	1.73-2.66
Control 2	821105	102	85.1-128
Panel 1	MF-101617-01	1.39	0.691-1.67
Panel 2	MF-101617-02	30.8	22.4-45.4
Panel 3	MF-101617-03	136	91.4-207

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	821105	5.62	3.54-7.91
Control 2	821105	483	409-614
Panel 1	MF-101617-01	4.69	3.06-6.42
Panel 2	MF-101617-02	32.1	17.1-50.9
Panel 3	MF-101617-03	366	242-551

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	821105	21.1	17.6-26.7
Control 2	821105	1105	906-1359
Panel 1	MF-101617-01	50.2	40.5-60.7
Panel 2	MF-101617-02	186	108-231
Panel 3	MF-101617-03	2675	2181-3314

UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	821105	131	90.6-138
Control 2	821105	11160	9059-13589
Panel 1	MF-101617-01	59.8	44.5-84.7
Panel 2	MF-101617-02	359	308-476
Panel 3	MF-101617-03	4951	4225-6337

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

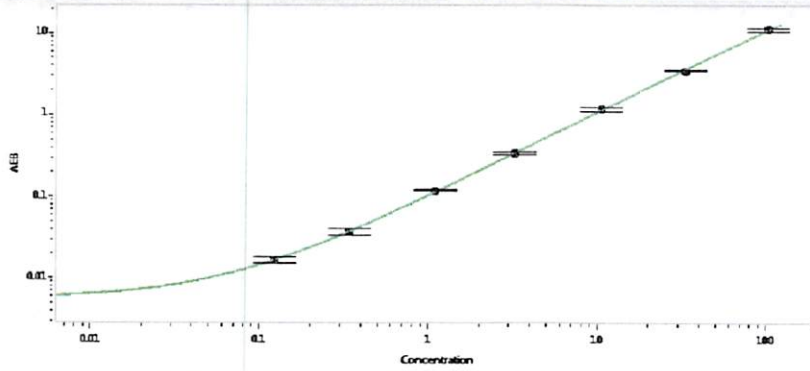
Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

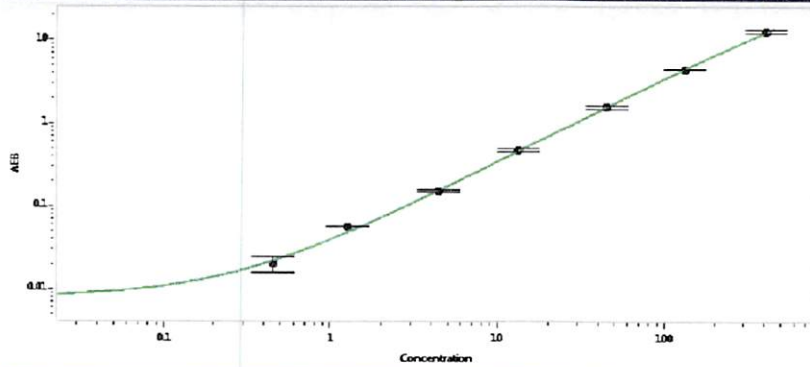
[Signature] 06 AUG 2018
Signature/Date

Tau Calibration Curve



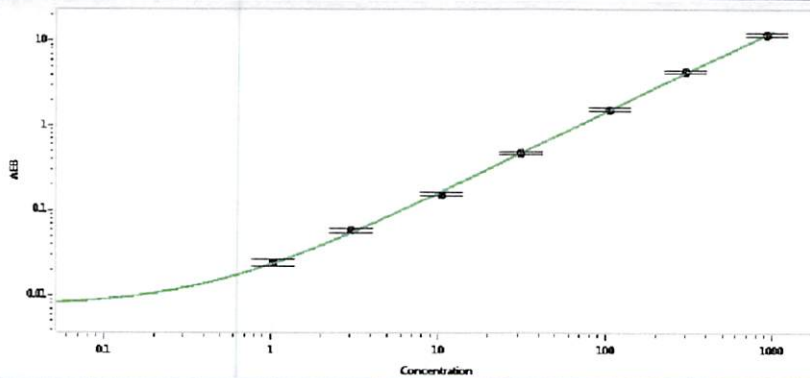
Calibrator Levels (pg/mL)	
A	0.000
B	0.124
C	0.341
D	1.10
E	3.24
F	10.5
G	33.0
H	103

NF-light® Calibration Curve



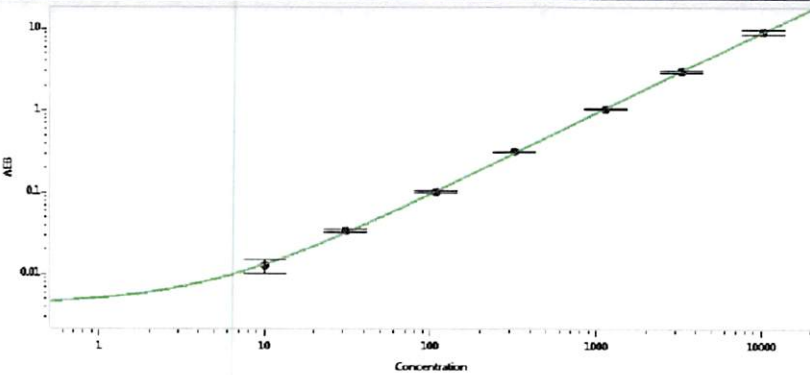
Calibrator Levels (pg/mL)	
A	0.000
B	0.451
C	1.27
D	4.39
E	13.1
F	44.6
G	132
H	406

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