

Product Number:	102153
Lot Number:	502101
Expiration:	28-Sep-2020
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102141	928015
Detector Reagent	102143	928003
SBG Reagent	102145	927307
Sample Diluent	102147	927105
Calibrators	102150	927601
Controls	102151-102152	927602
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is 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	927602	2.04	1.41-2.49
Control 2	927602	86.2	75.1-113
Panel 1	921101	0.630	0.309-2.15
Panel 2	921102	0.497	0.218-1.76
Panel 3	921103	19.7	13.7-40.3
Panel 4	921104	29.0	25.4-54.2

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	927602	6.48	5.34-8.01
Control 2	927602	507	445-668
Panel 1	921101	9.99	8.50-12.7
Panel 2	921102	8.68	6.95-10.4
Panel 3	921103	181	151-227
Panel 4	921104	630	560-840

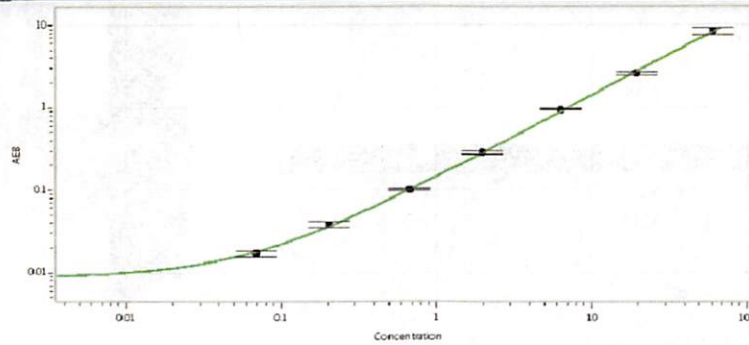
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	927602	23.0	18.3-27.5
Control 2	927602	1097	961-1441
Panel 1	921101	84.7	66.8-100
Panel 2	921102	60.9	45.9-68.9
Panel 3	921103	1177	899-1349
Panel 4	921104	6461	5080-8293

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	927602	76.6	54.0-115
Control 2	927602	10340	8035-12052
Panel 1	921101	82.6	72.7-139
Panel 2	921102	85.4	54.4-165
Panel 3	921103	2881	2680-4086
Panel 4	921104	7752	7104-11478

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

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	08 Nov 2019
	Name	Title	Signature/Date	

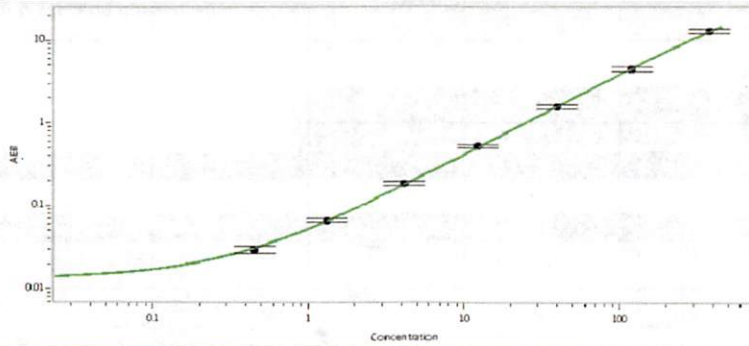
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.111
C	0.319
D	0.985
E	2.76
F	8.83
G	25.6
H	78.2

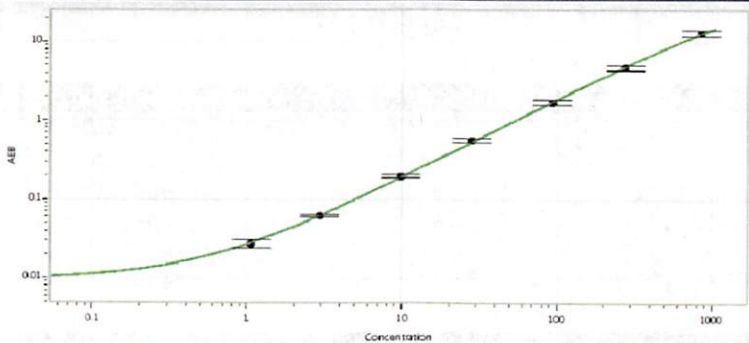
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.557
C	1.49
D	4.75
E	13.7
F	43.4
G	129
H	391

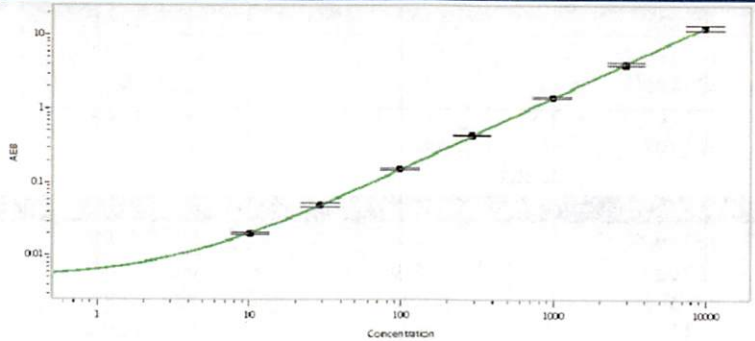
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.16
C	3.16
D	9.94
E	28.7
F	94.1
G	282
H	848

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	6.30
C	17.7
D	66.5
E	202
F	668
G	2032
H	6163