

Product Number:	102153
Lot Number:	502148
Expiration:	24-Jun-2020
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102141	931703
Detector Reagent	102143	931705
SBG Reagent	102145	917502
Sample Diluent	102147	930304
Calibrators	102150	931601 / 001304
Controls	102151-102152	931602
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	931602	2.11	1.67-2.51
Control 2	931602	96.2	77.8-117
Panel 1	921101	1.38	0.419-2.10
Panel 2	921102	1.14	0.311-1.73
Panel 3	921103	22.8	13.5-39.2
Panel 4	921104	39.6	26.8-52.8

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	6.85	5.44-8.16
Control 2	931602	627	491-737
Panel 1	921101	10.1	8.43-12.6
Panel 2	921102	9.10	7.00-10.5
Panel 3	921103	198	152-228
Panel 4	921104	734	565-847

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	22.3	17.1-25.7
Control 2	931602	1172	914-1371
Panel 1	921101	77.8	66.1-99.2
Panel 2	921102	55.6	45.7-68.5
Panel 3	921103	1323	924-1385
Panel 4	921104	7382	5201-8417

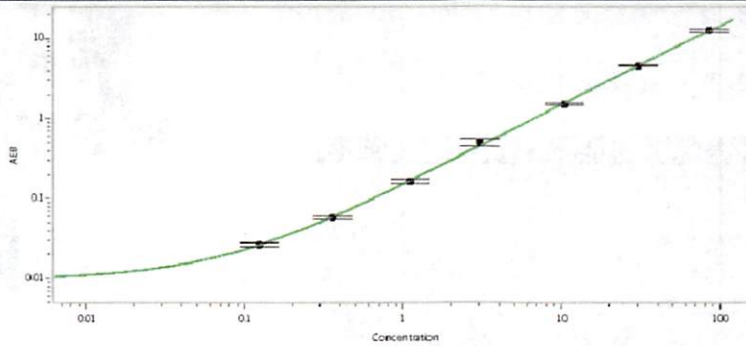
UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	145	118-187
Control 2	931602	14978	12112-18168
Panel 1	921101	82.5	64.3-140
Panel 2	921102	96.3	52.9-161
Panel 3	921103	3005	2589-4060
Panel 4	921104	8218	6857-11346

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

Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stodler 22 Jan 2020
Name Title Signature/Date

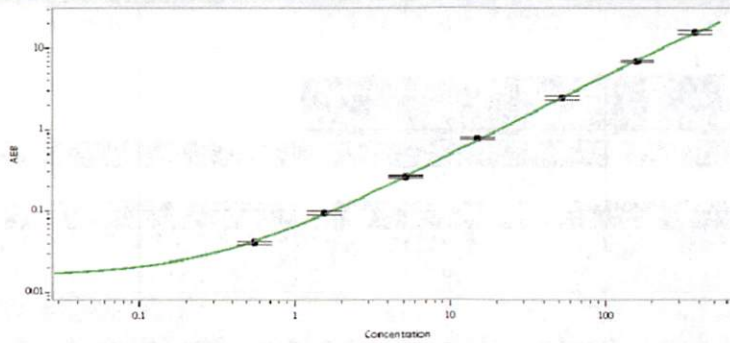
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.123
C	0.357
D	1.11
E	3.04
F	10.4
G	30.5
H	85.7

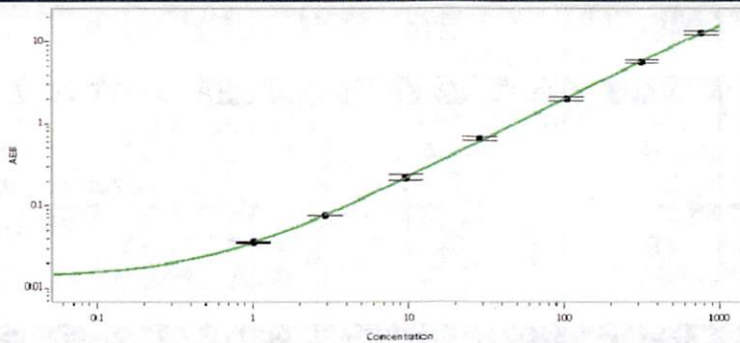
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.554
C	1.55
D	5.22
E	15.0
F	52.6
G	159
H	378

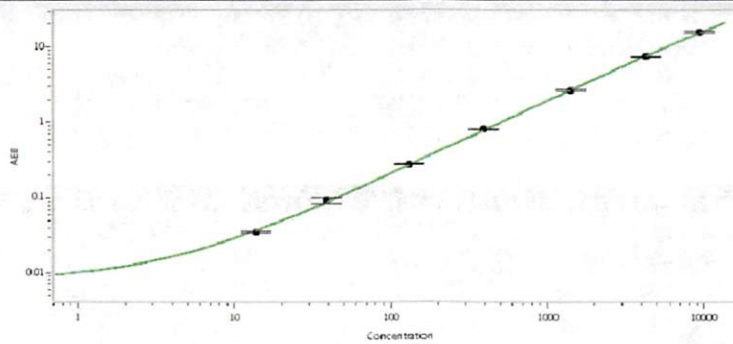
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.02
C	2.92
D	9.71
E	28.6
F	104
G	318
H	760

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	13.8
C	38.8
D	131
E	390
F	1413
G	4256
H	9366