

Product Number:	102153
Lot Number:	502219
Expiration:	30-Oct-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	003006
Sample Diluent	102147	930304
Calibrators	102150	931601 / 001304
Control 1	102151	931602
Control 2	102152	931602
Control 2	102152	934002
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ These reagents are not Kit Lot Specific

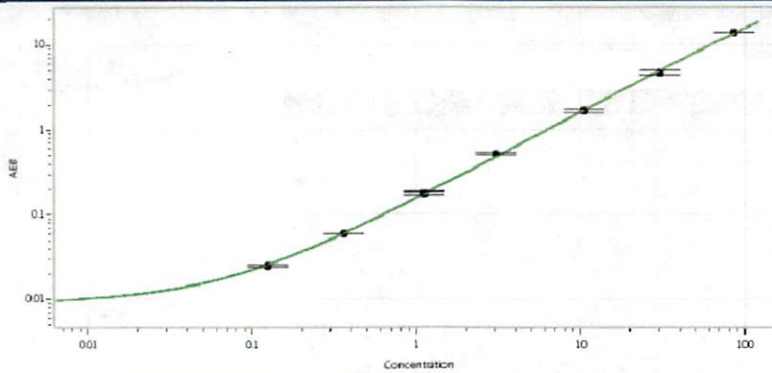
† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).

Release Materials			
Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	2.14	1.67-2.51
Control 2	931602	94.4	77.8-117
Control 2	934002	60.7	55.2-97.8
Panel 1	921101	1.50	0.419-2.10
Panel 2	921102	1.28	0.311-1.73
Panel 3	921103	22.6	13.5-39.2
Panel 4	921104	36.9	26.8-52.8
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	6.87	5.44-8.16
Control 2	931602	605	491-737
Control 2	934002	533	370-761
Panel 1	921101	10.0	8.43-12.6
Panel 2	921102	9.00	7.00-10.5
Panel 3	921103	193	152-228
Panel 4	921104	716	565-847
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	21.7	17.1-25.7
Control 2	931602	1126	914-1371
Control 2	934002	1054	779-1168
Panel 1	921101	78.0	66.1-99.2
Panel 2	921102	54.8	45.7-68.5
Panel 3	921103	1237	924-1385
Panel 4	921104	8093	5201-8417
UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931602	150	118-187
Control 2	931602	14106	12112-18168
Control 2	934002	12107	10243-15364
Panel 1	921101	88.1	64.3-140
Panel 2	921102	103	52.9-161
Panel 3	921103	2991	2589-4060
Panel 4	921104	7905	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
 Name: Patrizia Stadler Title: Assoc. Director, QC Signature/Date: P Stadler 08 Jul 2020

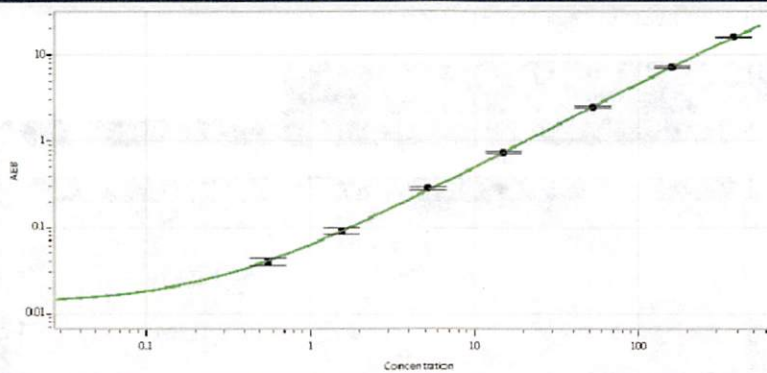
Tau Calibration Curve



**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

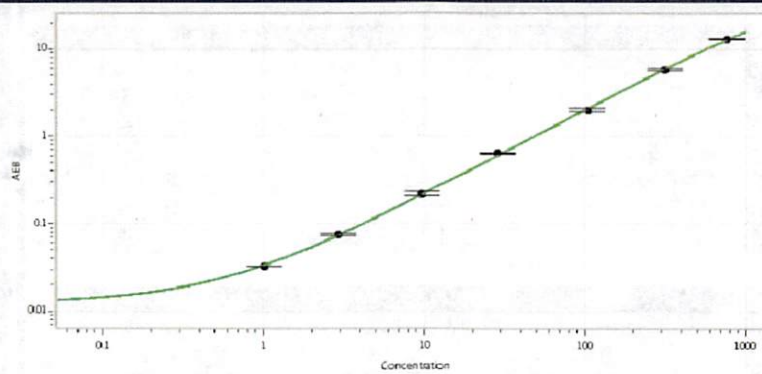
NF-light® Calibration Curve



**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

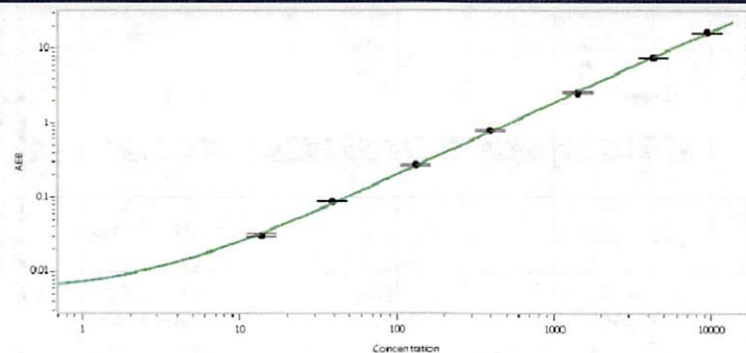
GFAP® Calibration Curve



**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