

Product Number:	103345
Lot Number:	503327
Expiration:	29-Dec-2022
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103338	210604
Detector Reagent	103339	212206
SBG Reagent	103340	136302
Sample Diluent	103341	201208
Calibrators	103342	134601
Control 1	103343	134602
Control 2	103344	134602
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex B SR-X Kit Instructions (KI-0123).

**Release Materials**

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	2.33	1.79-2.69
Control 2	134602	84.7	59.3-103
Panel 1	206807	12.6	10.8-16.1
Panel 2	206807	7.96	6.66-9.99
Panel 3	206807	26.0	21.3-32.0
Panel 4	206807	95.1	75.3-113

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	6.00	3.49-8.22
Control 2	134602	603	409-680
Panel 1	206807	9.75	7.64-11.5
Panel 2	206807	3.84	2.50-4.72
Panel 3	206807	153	98.1-164
Panel 4	206807	824	556-880

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	138	95.9-154
Control 2	134602	11595	8277-13641
Panel 1	206807	69.8	50.2-82.4
Panel 2	206807	27.3	12.4-33.2
Panel 3	206807	977	762-1143
Panel 4	206807	5075	3908-5861

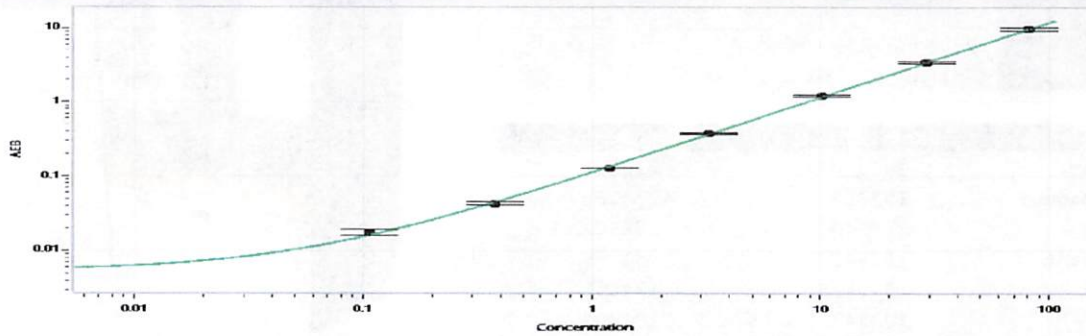
UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	110	67.4-150
Control 2	134602	8340	6654-9981
Panel 1	206807	65.0	42.6-88.4
Panel 2	206807	62.6	37.9-82.4
Panel 3	206807	962	790-1185
Panel 4	206807	6604	5493-9112

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

Review/Approval Sonia Khilkin QC Manager Sonia Khilkin / 13 JUL 2022  
 Name Title Signature/Date



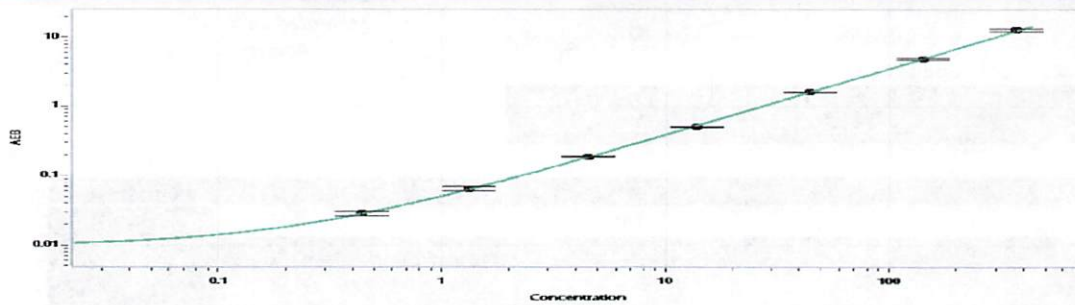
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.107
C	0.379
D	1.19
E	3.22
F	10.07
G	29.2
H	80.6

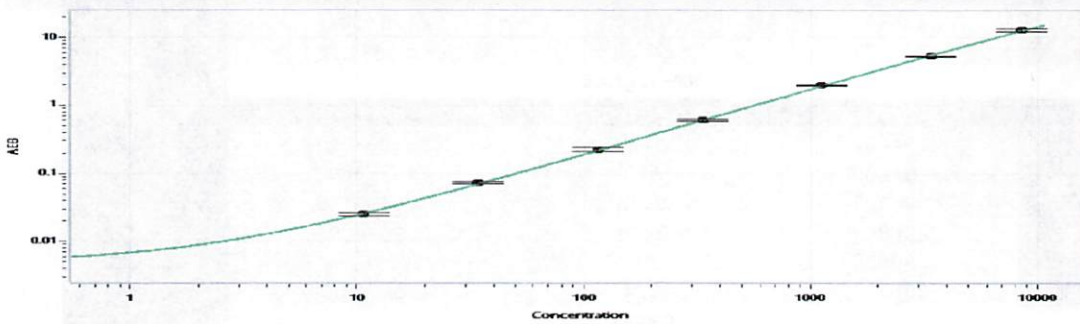
**NF-light® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.441
C	1.33
D	4.53
E	13.8
F	44.4
G	142
H	372

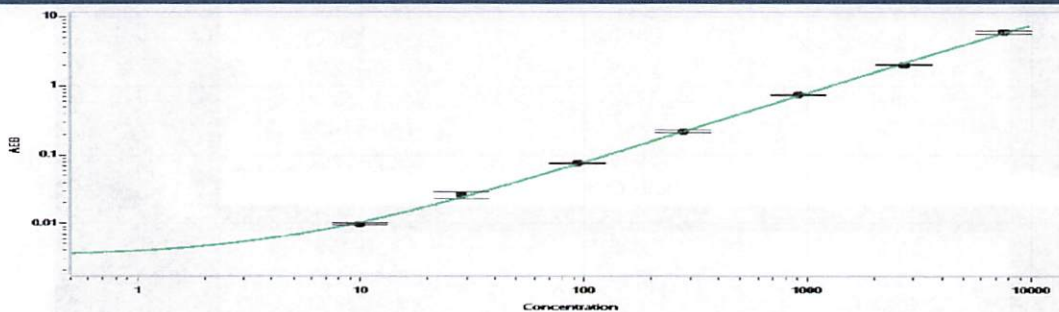
**GFAP® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	10.77
C	34.2
D	115.7
E	334
F	1129
G	3385
H	8525

**UCHL1® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	9.9
C	28.2
D	93.7
E	278
F	909
G	2720
H	7525