

Product Number:	103345
Lot Number:	502123
Expiration:	18-Nov-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	935008
Detector Reagent	103339	001413
SBG Reagent	103340	932204
Sample Diluent	103341	932203
Calibrators	103342	931901
Controls	103343-103344	931902
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).



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

**Release Materials**

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	1.96	1.42-2.30
Control 2	931902	84.9	70.4-106
Panel 2	921102	0.898	0.557-2.95
Panel 3	921103	31.1	22.4-33.6
Panel 4	921104	47.9	36.4-58.8

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	6.06	4.83-7.25
Control 2	931902	530	440-661
Panel 2	921102	7.18	5.74-8.61
Panel 3	921103	201	145-217
Panel 4	921104	663	542-813

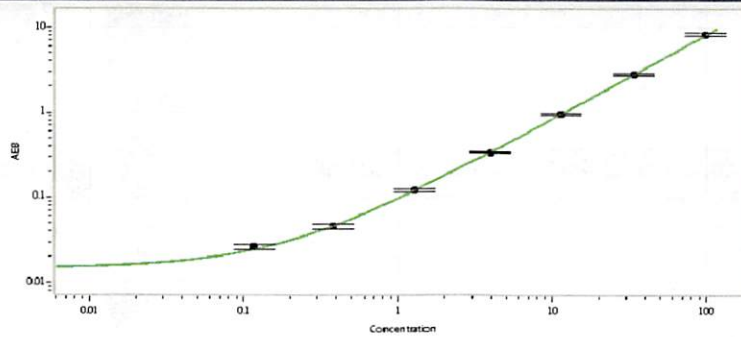
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	91.3	73.0-109
Control 2	931902	9876	8070-12105
Panel 2	921102	43.3	27.5-53.5
Panel 3	921103	1128	625-1222
Panel 4	921104	6198	3418-6873

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	131	87.8-132
Control 2	931902	11830	8175-12262
Panel 2	921102	112	84.3-127
Panel 3	921103	3924	2320-4225
Panel 4	921104	9768	6239-10760

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>	<i>22 Jan 2020</i>
	Name	Title	Signature/Date	

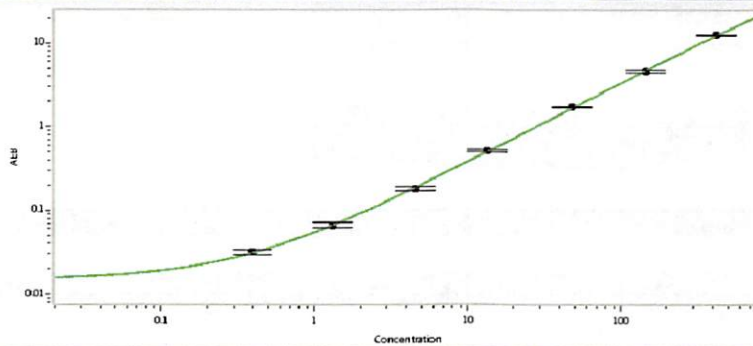
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.117
C	0.377
D	1.28
E	3.95
F	11.4
G	33.7
H	97.8

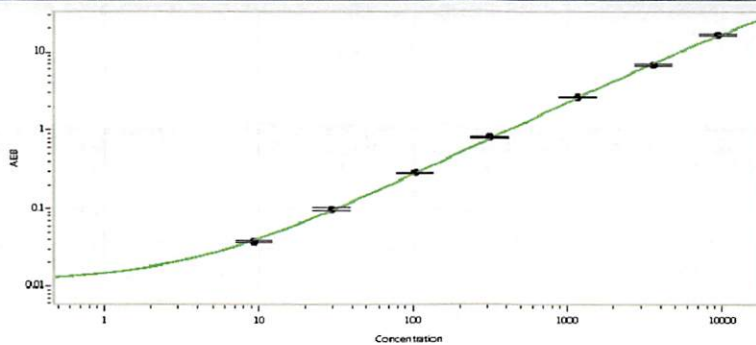
**NF-light® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.393
C	1.32
D	4.58
E	13.3
F	47.9
G	146
H	421

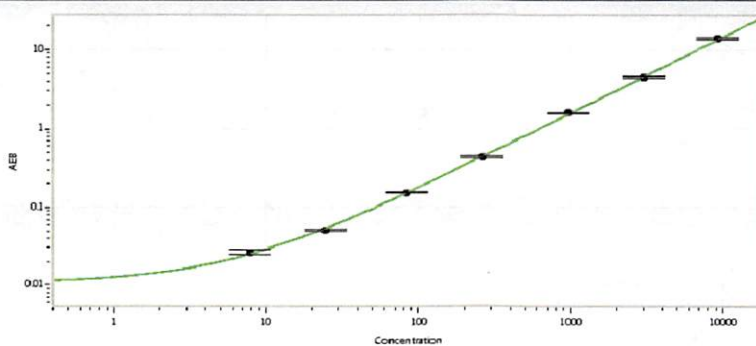
**GFAP® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	9.33
C	29.4
D	104
E	314
F	1165
G	3596
H	9411

**UCHL1® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	7.89
C	24.4
D	83.8
E	260
F	971
G	3007
H	9266