

Product Number:	103345
Lot Number:	501915
Expiration:	19-Feb-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	919211
Detector Reagent	103339	922103
SBG Reagent	103340	905403
Sample Diluent	103341	905010
Calibrators	103342	921401
Controls	103343-103344	921402
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>



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

<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	921401	1.60	1.25-1.87
Control 2	921402	78.2	69.4-104
Panel 1	921101	2.81	1.93-3.29
Panel 2	921102	2.25	1.77-2.68
Panel 3	921103	28.3	22.8-34.2
Panel 4	921104	47.6	39.3-58.9

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921401	5.60	4.52-6.78
Control 2	921402	567	451-677
Panel 1	921101	8.25	7.11-10.7
Panel 2	921102	7.19	5.62-8.42
Panel 3	921103	189	138-208
Panel 4	921104	744	550-824

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921401	97.5	78.5-118
Control 2	921402	10299	8524-12786
Panel 1	921101	53.6	43.9-68.1
Panel 2	921102	41.7	30.4-47.3
Panel 3	921103	908	657-985
Panel 4	921104	5056	3715-5572

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921401	116	65.1-134
Control 2	921402	10126	7481-11322
Panel 1	921101	92.0	64.9-120
Panel 2	921102	111	80.1-122
Panel 3	921103	3018	2391-3587
Panel 4	921104	7856	6249-9374

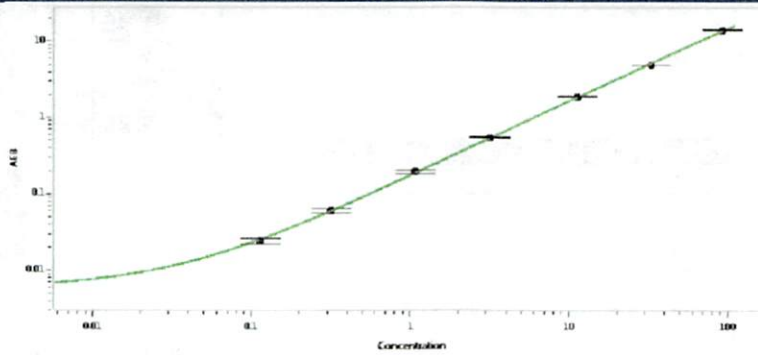
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval: Patrizia Stadler Assoc. Director, QC  
Name: Patrizia Stadler Title: Assoc. Director, QC

P. Stadler 12 Aug 2019  
Signature/Date



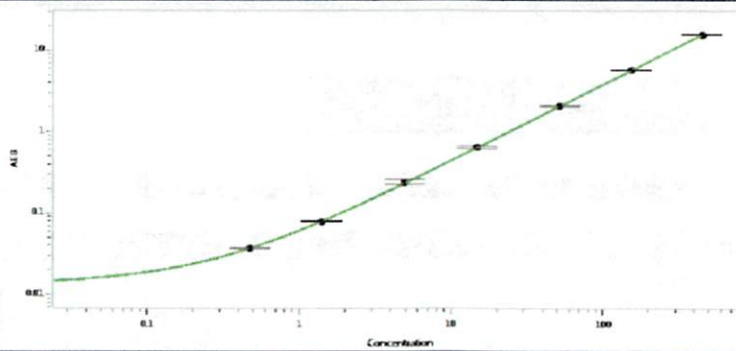
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.113
C	0.317
D	1.07
E	3.17
F	11.3
G	32.9
H	92.8

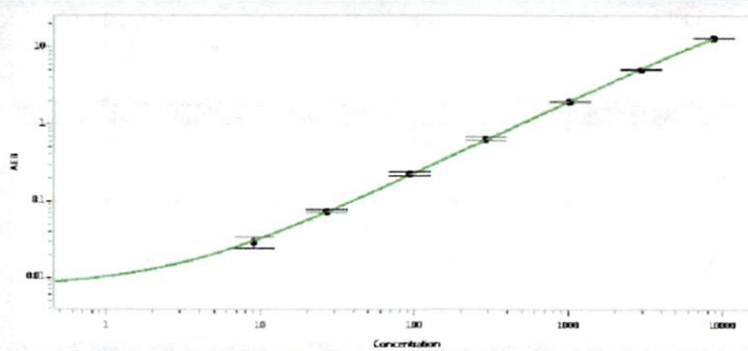
**NF-light® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.479
C	1.41
D	4.92
E	14.7
F	52.1
G	153
H	448

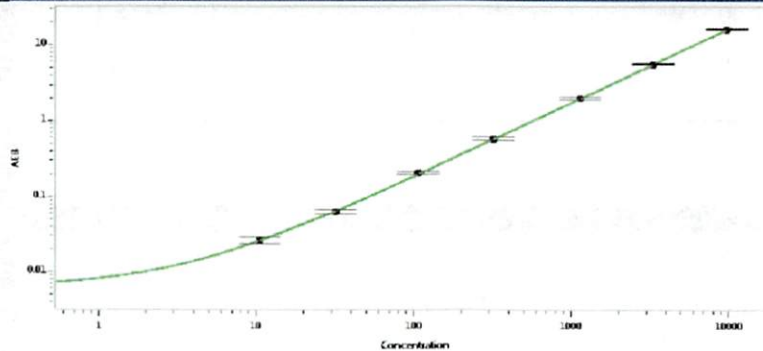
**GFAP® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	9.08
C	26.7
D	93.1
E	288
F	1012
G	2976
H	8671

**UCHL1® Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	10.6
C	32.1
D	108
E	322
F	1146
G	3332
H	9833