

Product Number:	102153
Lot Number:	502600
Expiration:	24-Aug-2021
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102141	030708
Detector Reagent	102143	031506
SBG Reagent	102145	028313
Sample Diluent	102147	030004
Calibrators	102150	019901
Control 1	102151	019902
Control 2	102152	019902
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 A SR-X Kit Instructions (KI-0031).



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

**Release Materials**

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	2.43	1.62-2.68
Control 2	019902	85.7	69.7-105
Panel 1	921101	1.37	0.419-2.10
Panel 2	921102	1.10	0.311-1.73
Panel 3	921103	25.4	14.4-38.7

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	6.72	5.18-7.77
Control 2	019902	535	428-642
Panel 1	921101	9.01	8.43-12.6
Panel 2	921102	7.60	7.00-10.5
Panel 3	921103	200	151-240

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	29.8	14.1-33.1
Control 2	019902	989	824-1235
Panel 1	921101	76.7	66.1-99.2
Panel 2	921102	53.3	45.7-68.5
Panel 3	921103	1104	927-1391

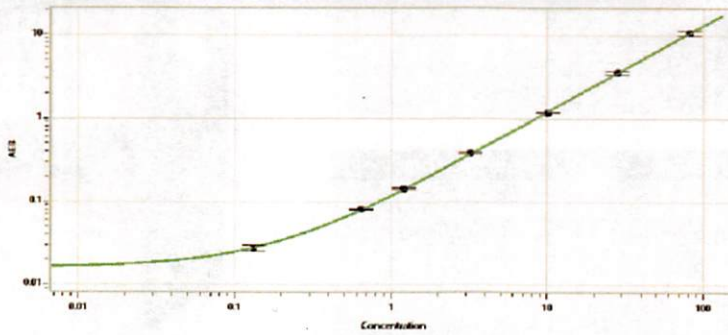
UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	117	81.8-132
Control 2	019902	8589	7118-10677
Panel 1	921101	106	64.3-140
Panel 2	921102	105	52.9-161
Panel 3	921103	3660	1662-5492

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 Title

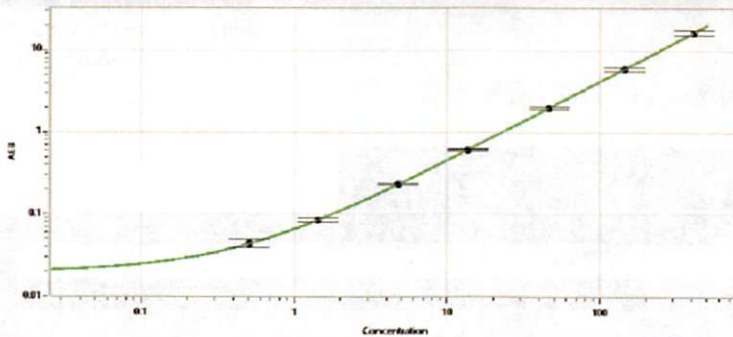
P. Stadler 18 Nov 2020  
Signature/Date

**Tau Calibration Curve**



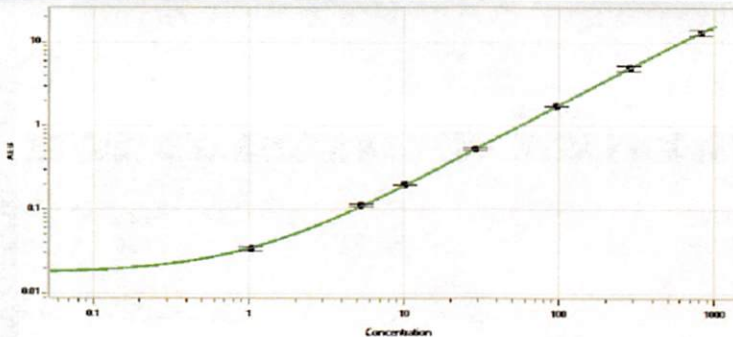
Calibrator Levels (pg/mL)	
A	0.000
B	0.133
C	0.647
D	1.22
E	3.22
F	10.1
G	27.9
H	81.1

**NF-light® Calibration Curve**



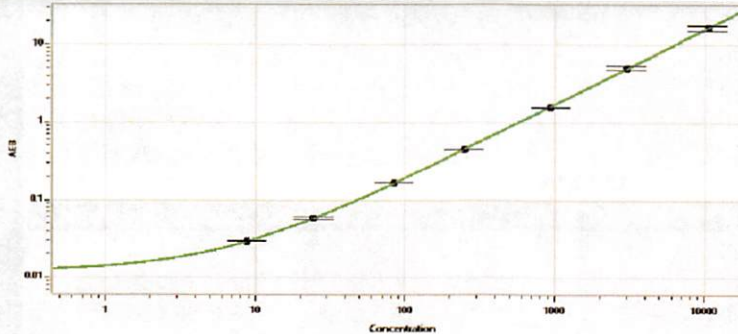
Calibrator Levels (pg/mL)	
A	0.000
B	0.506
C	1.43
D	4.81
E	13.7
F	46.4
G	144
H	410

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	1.04
C	5.29
D	10.1
E	28.8
F	97.0
G	282
H	820

**UCHL1® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	8.77
C	24.4
D	84.4
E	248
F	940
G	2998
H	10508