

Product Number:	103345
Lot Number:	502538
Expiration:	4-Mar-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	006410
Detector Reagent	103339	006413
SBG Reagent	103340	006502
Sample Diluent	103341	006503
Calibrators	103342	921401
Control 1	103343	921402
Control 2	103344	921402
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	921402	1.44	1.24 - 1.96
Control 2	921402	77.4	67.0 - 101
Panel 2	921102	0.999	0.312 - 2.93
Panel 3	921103	23.0	20.3 - 34.0
Panel 4	921104	35.2	29.9 - 61.1

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921402	6.19	4.35 - 7.64
Control 2	921402	606	457 - 686
Panel 2	921102	7.26	5.75 - 8.63
Panel 3	921103	199	147 - 221
Panel 4	921104	775	555 - 832

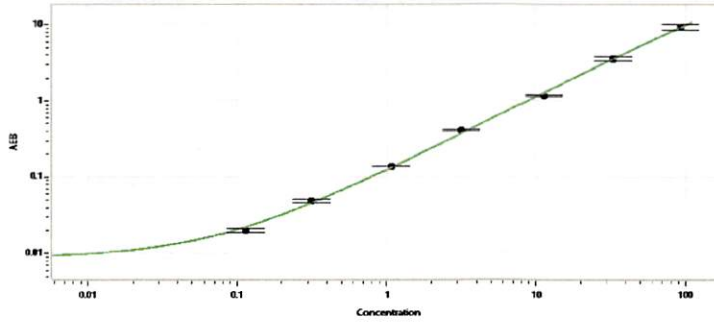
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921402	94.8	77.9 - 117
Control 2	921402	10909	8606 - 12908
Panel 2	921102	40.5	27.8 - 53.1
Panel 3	921103	883	638 - 1195
Panel 4	921104	5394	3584 - 6789

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921402	128	68.0 - 146
Control 2	921402	11776	7227 - 13217
Panel 2	921102	106	84.5 - 127
Panel 3	921103	3297	2398 - 4156
Panel 4	921104	8992	6436 - 10728

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

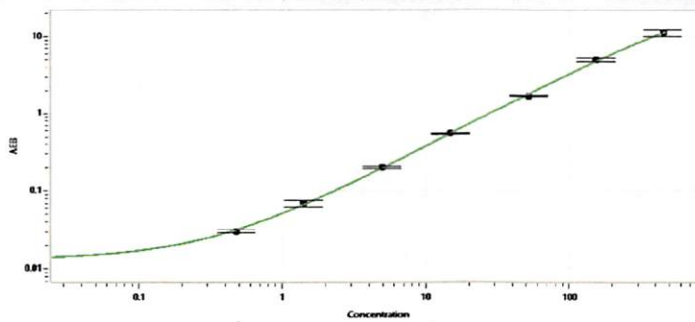
Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	30 Sep 2020
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	

**Tau Calibration Curve**



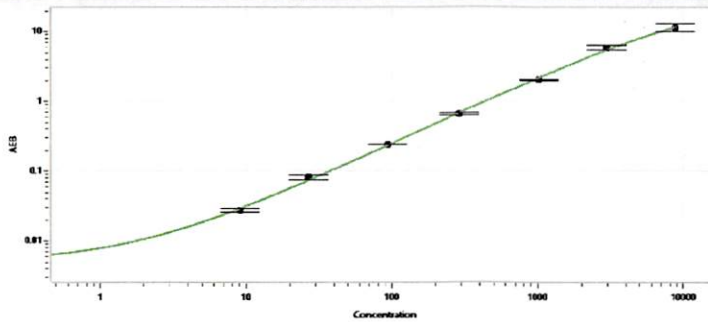
Calibrator Levels (pg/mL)	
A	0.000
B	0.113
C	0.317
D	1.07
E	3.17
F	11.3
G	32.9
H	92.8

**NF-light® Calibration Curve**



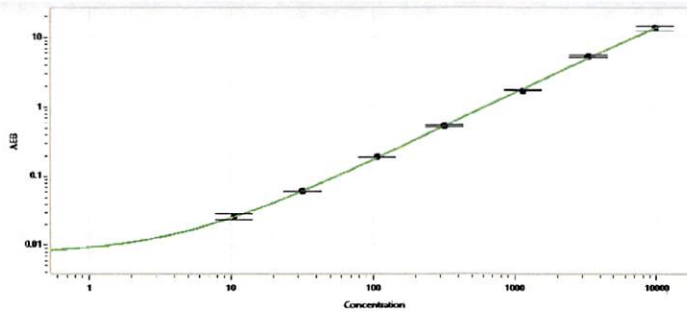
Calibrator Levels (pg/mL)	
A	0.000
B	0.479
C	1.41
D	4.92
E	14.7
F	52.1
G	153
H	448

**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