

Product Number:	103345
Lot Number:	503270
Expiration:	10-Nov-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	201207
Detector Reagent	103339	211805
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 ¹
RGP Reagent	103159	N/A ¹

¹ 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	134602	2.19	1.77-2.68
Control 2	134602	85.2	56.6-104
Panel 1	206807	14.2	10.9-16.4
Panel 2	206807	8.95	6.73-10.1
Panel 3	206807	27.1	21.4-32.1
Panel 4	206807	92.6	75.1-113

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	6.06	3.25-8.39
Control 2	134602	546	407-653
Panel 1	206807	9.62	7.60-11.4
Panel 2	206807	3.72	2.55-4.56
Panel 3	206807	118	100-151
Panel 4	206807	655	553-830

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	115	96.7-146
Control 2	134602	11034	8047-13553
Panel 1	206807	70.3	48.6-82.3
Panel 2	206807	23.1	13.8-29.5
Panel 3	206807	958	757-1136
Panel 4	206807	4752	3870-5804

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134602	101	63.2-154
Control 2	134602	8051	6650-9975
Panel 1	206807	67.6	41.8-89.5
Panel 2	206807	55.7	39.3-79.7
Panel 3	206807	951	794-1192
Panel 4	206807	6479	5730-9224

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

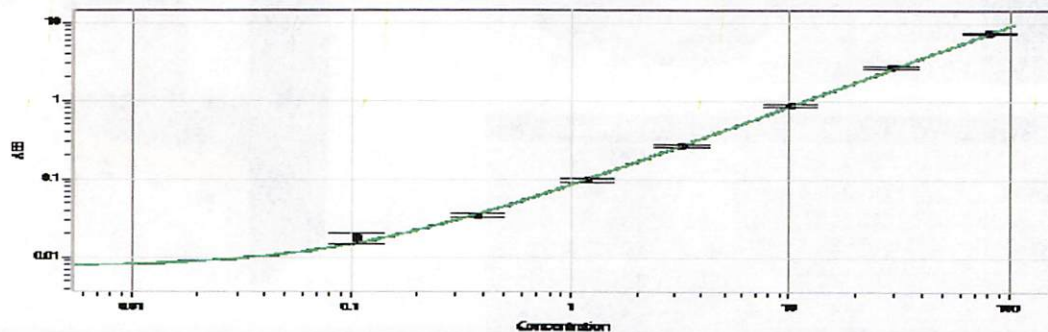
Review/Approval

Matt Smart
Name

QC Director
Title

M. A. 06 Jun 2022
Signature/Date

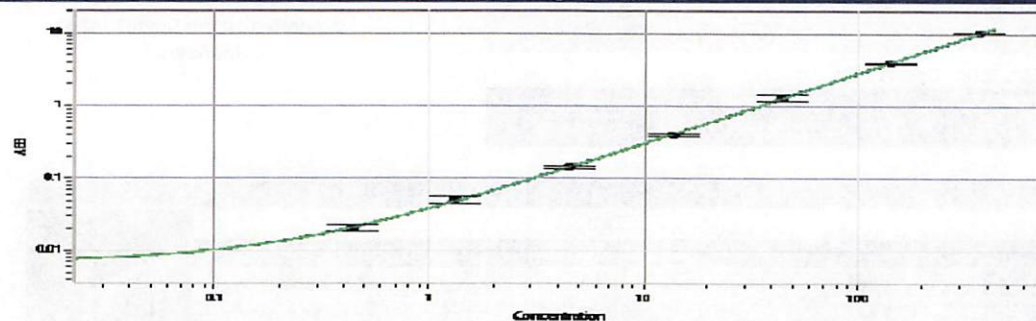
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.107
C	0.379
D	1.19
E	3.22
F	10.1
G	29.2
H	80.6

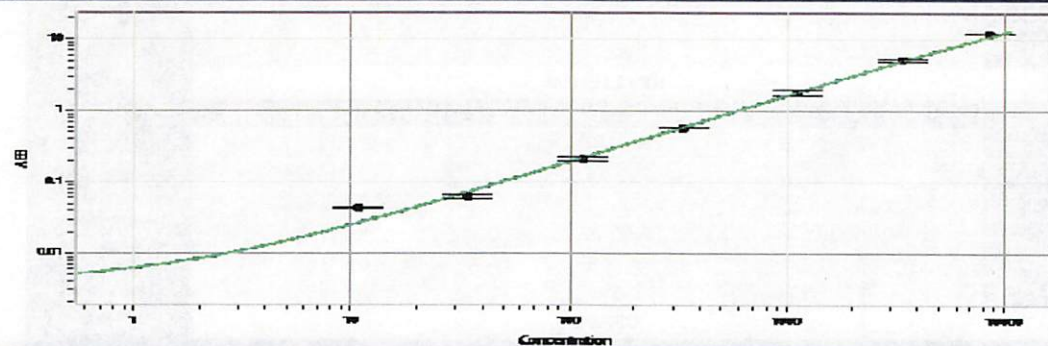
NF-light® Calibration Curve



**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

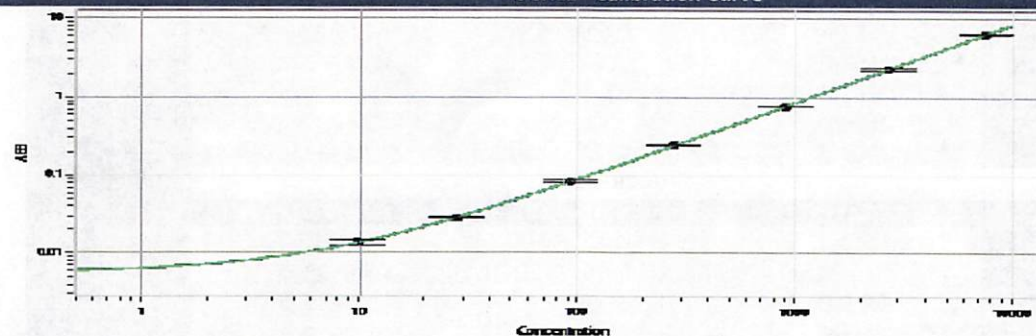
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	10.8
C	34.2
D	116
E	334
F	1129
G	3385
H	8525

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	9.93
C	28.2
D	93.7
E	278
F	909
G	2720
H	7525