

Product Number:	102153
Lot Number:	502410
Expiration:	21-Jan-2021
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102141	015001
Detector Reagent	102143	015002
SBG Reagent	102145	003006
Sample Diluent	102147	002117
Calibrators	102150	015301
Control 1	102151	015302
Control 2	102152	015302
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 A SR-X Kit Instructions (KI-0031).

Release Materials

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	015302	2.38	1.74-2.74
Control 2	015302	90.8	74.4-111
Panel 2	921102	0.490	0.311-1.73
Panel 3	921103	20.2	14.4-38.7
Panel 4	921104	30.9	27.7-53.0

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	015302	6.82	5.36-8.04
Control 2	015302	531	433-649
Panel 2	921102	7.73	7.00-10.5
Panel 3	921103	174	151-240
Panel 4	921104	620	561-897

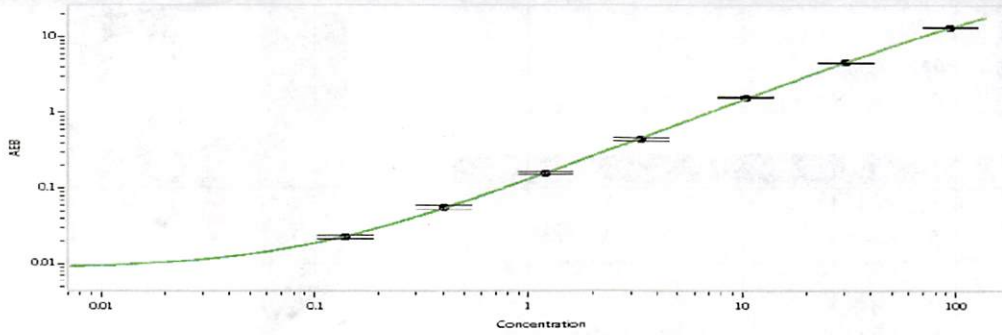
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	015302	22.8	17.8-26.7
Control 2	015302	1106	888-1332
Panel 2	921102	56.7	45.7-68.5
Panel 3	921103	1097	927-1391
Panel 4	921104	7832	3970-8978

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	015302	108	75.4-124
Control 2	015302	9619	7503-11255
Panel 2	921102	127	52.9-161
Panel 3	921103	3581	1662-5492
Panel 4	921104	9703	4277-15688

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

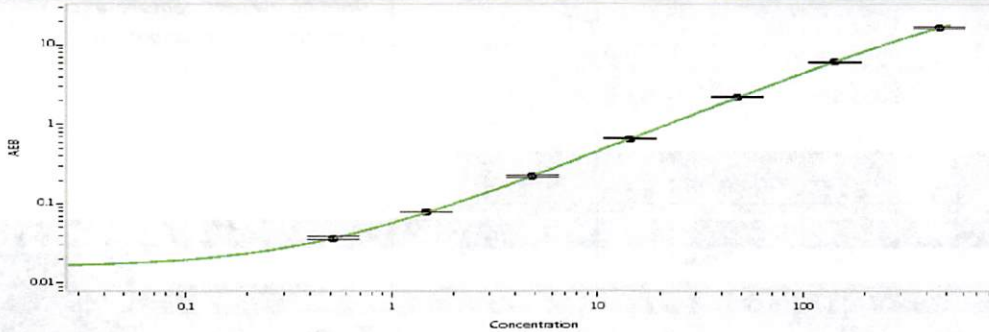
Review/Approval Shivani Goel QC Manager Shivani Goel 15 Jul 2020
Name Title Signature/Date

Tau Calibration Curve



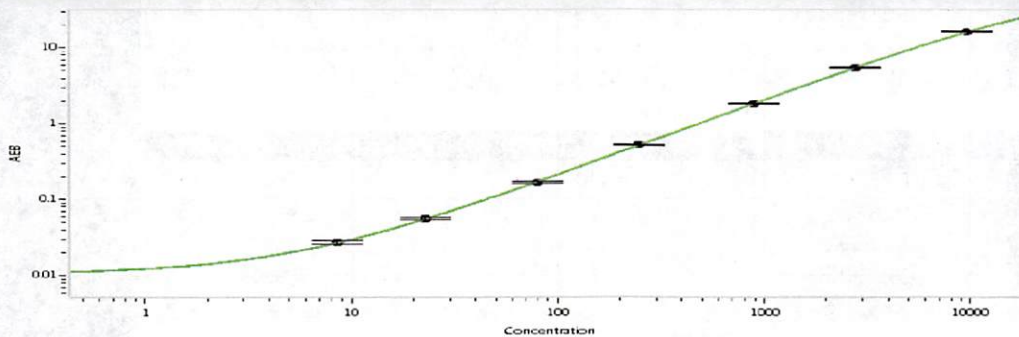
Calibrator Levels (pg/mL)	
A	0.000
B	0.139
C	0.400
D	1.20
E	3.34
F	10.4
G	30.6
H	94.1

NF-light® Calibration Curve



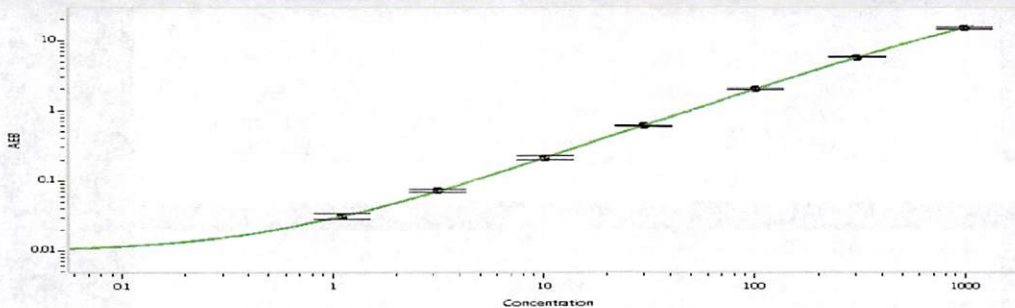
Calibrator Levels (pg/mL)	
A	0.000
B	0.522
C	1.48
D	4.83
E	14.4
F	48.1
G	143
H	461

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.11
C	3.17
D	10.2
E	29.8
F	101
G	307
H	991

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	8.51
C	22.9
D	79.5
E	246
F	886
G	2769
H	9667