

Neuro 2-Plex B Bulk Kit	
Product Number:	104080
Lot Number:	503398
Expiration:	15-Feb-2023
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	204906
Detector Reagent	103512	204907
SBG Reagent	103514	205001
Sample Diluent	103516	205002
Calibrators	103517	135301
Control 0	104079	205809
Control 1	104078	205809
Control 2	104077	205809
RGP Reagent ¹	103159	N/A

¹ RGP is not Kit Lot Specific

Release Materials			
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 0	205809	23.7	18.4 - 27.5
Control 1	205809	92.9	72.3 - 108
Control 2	205809	283	223 - 334
Panel 1	207003	13.7	10.6 - 15.9
Panel 2	207003	6.89	4.58 - 7.60
Panel 3	207003	130	101 - 151
Panel 4	207003	379	284 - 427
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 0	205809	89.5	30.0 - 103
Control 1	205809	505	204 - 569
Control 2	205809	2499	1088 - 2840
Panel 1	207003	67.3	55.3 - 83.7
Panel 2	207003	77.1	62.0 - 102
Panel 3	207003	902	749 - 1124
Panel 4	207003	5865	4803 - 7205

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

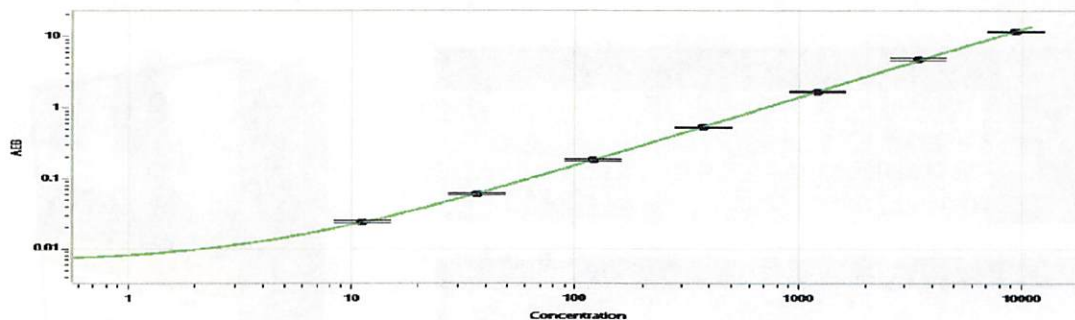
Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 22 Mar 2022
Signature/Date

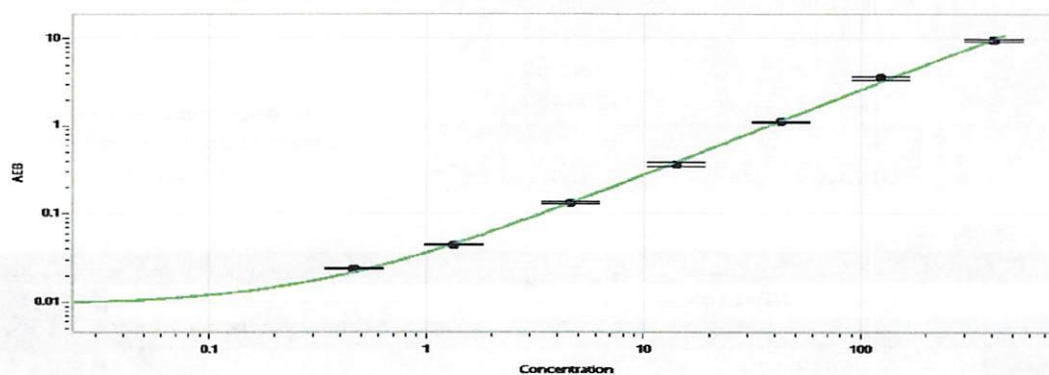
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	11.2
C	36.5
D	121
E	373
F	1224
G	3440
H	9416

NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.468
C	1.33
D	4.57
E	14.1
F	42.5
G	122
H	410