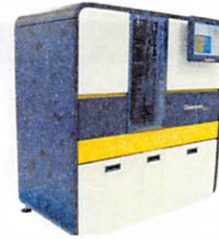


Neuro 2-Plex B Kit	
Product Number:	103520
Lot Number:	503343
Expiration:	4-Oct-2022
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	127706
Detector Reagent	103512	127802
SBG Reagent	103514	127703
Sample Diluent	103516	127704
Calibrators	103517	203302
Control 1	103518	203303
Control 2	103519	203303
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP / SBG Diluent (SR-X) are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa® Neuro 2-Plex B SR-X Kit Instructions (KI-0062).



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

Release Materials

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203303	5.65	4.53-7.10
Control 2	203303	538	435-653
Panel 1	207003	13.3	10.7-16.1
Panel 2	207003	6.5	4.77-7.68
Panel 3	207003	125	101-151
Panel 4	207003	324	273-416

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203303	92.8	72.2-111
Control 2	203303	9247	7332-10997
Panel 1	207003	64.7	49.4-83.9
Panel 2	207003	79.2	52.9-104
Panel 3	207003	859	684-1104
Panel 4	207003	5559	4259-7215

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

Review/Approval

Shivani Goel

QC Manager

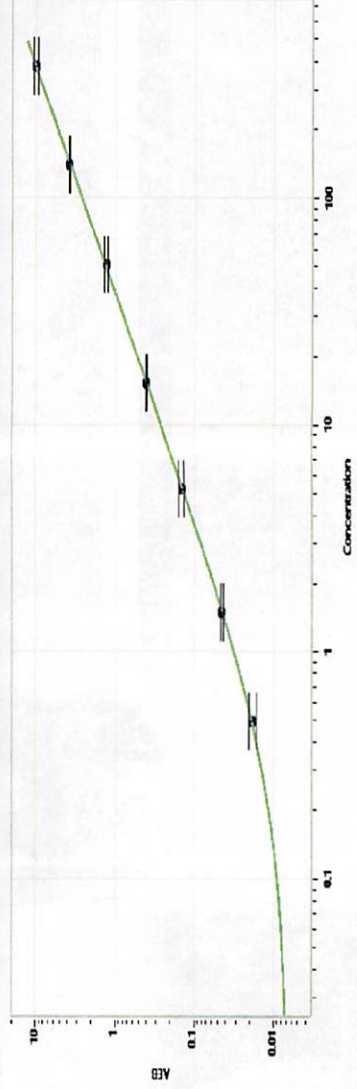
Shivani Goel 29 Mar 2022

Name

Title

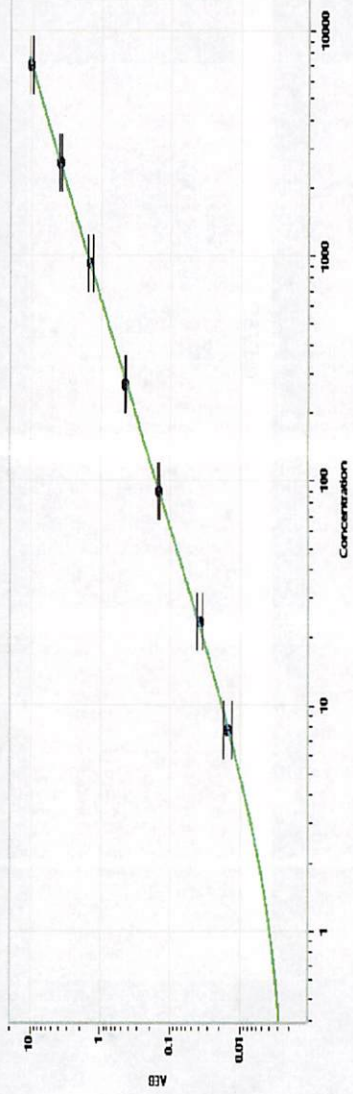
Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.492
C	1.50
D	5.20
E	15.3
F	50.7
G	140
H	377

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	7.75
C	23.6
D	89.5
E	268
F	925
G	2601
H	7092