

Neuro 4-Plex E Kit	
Product Number:	103670
Lot Number:	502932
Expiration:	14-May-2022
Neuro 4-Plex E Bulk Kit	
Product Number:	103971
Lot Number:	502959
Platform(s) :	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	113408
Detector Reagent	103655	113409
SBG Reagent	103657	113417
Plasma Sample Diluent	103659	113418
CSF Sample Diluent ¹	103727	113416
Calibrators	103667	114114
Control 1	103668	114115
Control 2	103669	114115
RGP Reagent ²	103159	N/A

¹ CSF Sample Diluent is included only in Neuro 4-Plex E kit (Product Number 103670)

² RGP is not Kit Lot Specific

Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114115	17.0	12.4-18.5
Control 2	114115	120	93.0-140
Panel 1	027301	29.7	23.3-34.9
Panel 2	018109	52.5	42.3-64.6
Panel 3	018109	134	108-162
Abeta 42			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114115	6.10	4.33-6.50
Control 2	114115	31.3	23.7-35.6
Panel 1	027301	2.25	2.00-3.00
Panel 2	018109	20.3	16.6-24.8
Panel 3	018109	42.6	36.0-54.0
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114115	179	142-212
Control 2	114115	3930	2966-4449
Panel 1	027301	45.2	37.5-56.2
Panel 2	018109	559	446-668
Panel 3	018109	2347	1916-2948
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114115	27.2	18.6-27.8
Control 2	114115	583	416-629
Panel 1	027301	10.7	7.10-12.6
Panel 2	018109	79.4	59.5-89.3
Panel 3	018109	432	310-474

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

Review/Approval

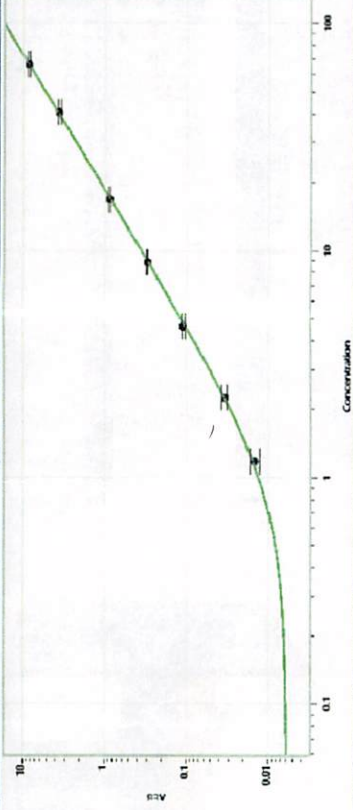
Richa Shrestha
Name

QC Supervisor
Title

Signature/Date

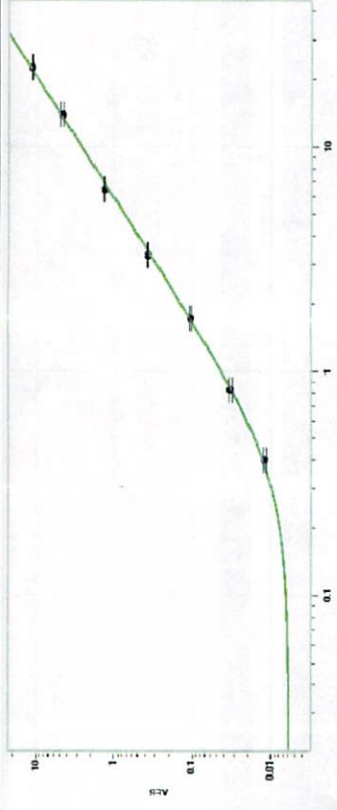
[Signature] 29 APR 2022

Abeta 40 Calibration Curve



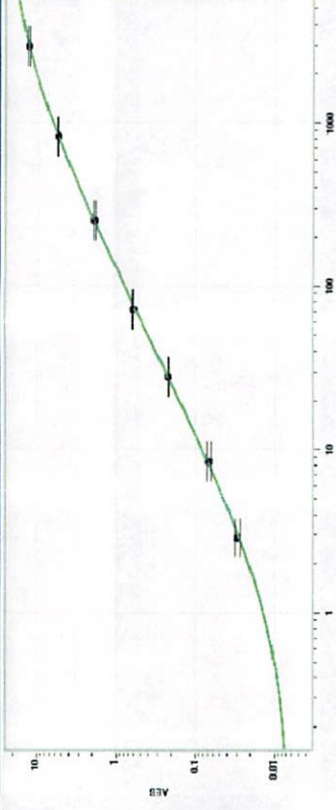
Calibrator Levels (pg/ml)	
A	0.000
B	1.18
C	2.25
D	4.63
E	8.79
F	16.7
G	40.4
H	65.5

Abeta 42 Calibration Curve



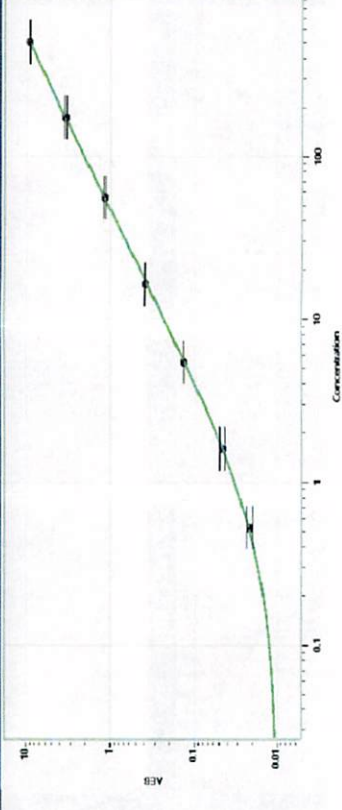
Calibrator Levels (pg/ml)	
A	0.000
B	0.403
C	0.825
D	1.72
E	3.29
F	6.43
G	13.8
H	22.4

GFAP® Calibration Curve



Calibrator Levels (pg/ml)	
A	0.000
B	5.01
C	15.3
D	50.8
E	139
F	479
G	1554
H	5204

NF-light® Calibration Curve



Calibrator Levels (pg/ml)	
A	0.000
B	0.525
C	1.60
D	5.47
E	16.2
F	55.3
G	173
H	495