

Product Number:	103670
Lot Number:	503105
Expiration:	10-Aug-2022
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	122801
Detector Reagent	103655	122802
SBG Reagent	103657	122815
Plasma Sample Diluent	103659	122813
CSF Sample Diluent	103727	122814
Calibrators	103667	123507
Control 1	103668	123508
Control 2	103669	123508
RGP Reagent	103159	N/A

¹ RGP is not Kit Lot Specific


Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	123508	12.5	8.68-17.0
Control 2	123508	87.3	63.2-120
Panel 1	027301	24.3	23.3 - 34.9
Panel 3	018109	115	100 - 163
Abeta 42			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	123508	4.83	4.15-6.22
Control 2	123508	29.9	24.3-36.4
Panel 1	027301	2.04	2.00 - 3.00
Panel 3	018109	43.8	36.0 - 54.0
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	123508	184	146-219
Control 2	123508	3639	3033-4549
Panel 1	027301	43.8	37.5 - 56.2
Panel 3	018109	2189	1916 - 2948
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	123508	17.3	15.1-22.7
Control 2	123508	395	333-500
Panel 1	027301	9.18	7.10 - 12.6
Panel 3	018109	412	310 - 474

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

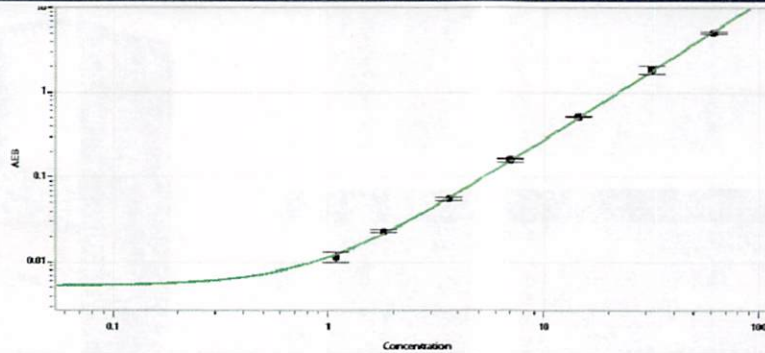
Review/Approval

Junpil Park
Name

QC Supervisor
Title

 07 Sep 2021
Signature/Date

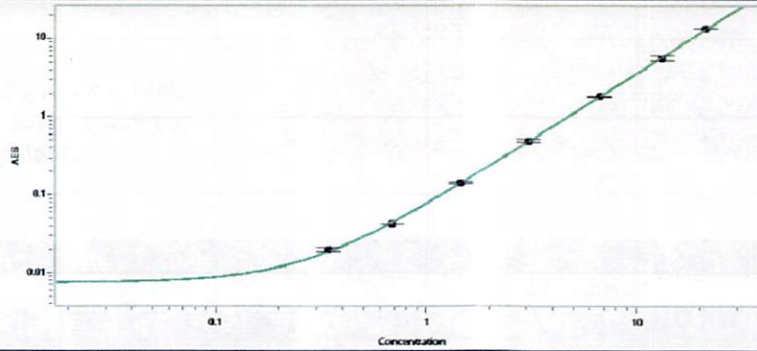
Abeta 40 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.08
C	1.80
D	3.62
E	7.00
F	14.6
G	31.7
H	61.5

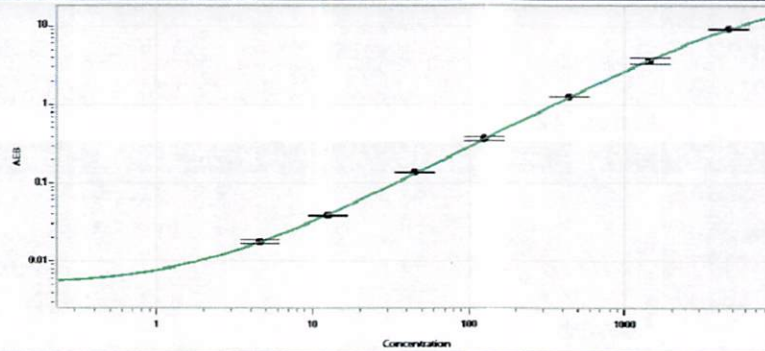
Abeta 42 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.343
C	0.690
D	1.46
E	3.05
F	6.63
G	13.1
H	21.1

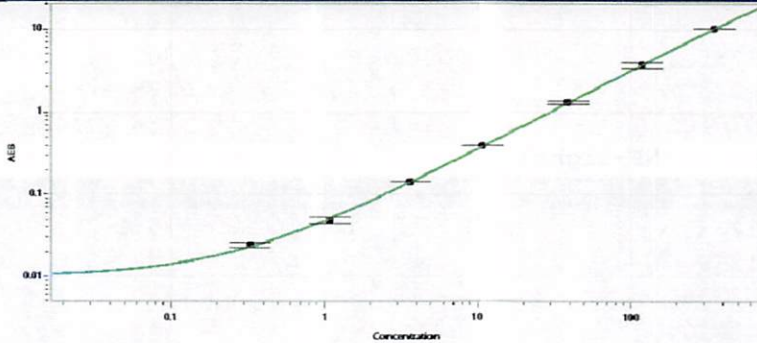
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	4.57
C	12.5
D	45.7
E	126
F	445
G	1443
H	4652

NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.330
C	1.09
D	3.64
E	10.7
F	39.1
G	118
H	352