

Product Number:	103670
Lot Number:	503237
Expiration:	18-Sep-2022
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	126101
Detector Reagent	103655	128801
SBG Reagent	103657	126105
Plasma Sample Diluent	103659	126202
CSF Sample Diluent	103727	126203
Calibrators	103667	132501
Control 1	103668	132502
Control 2	103669	132502
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	132502	22.9	12.9-25.0
Control 2	132502	139	105-157
Panel 1	027301	30.7	20.9-35.9
Panel 2	124507	97.0	69.2-104
Panel 3	125905	130	92.1-138
Abeta 42			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132502	6.57	4.73-7.10
Control 2	132502	36.2	26.7-40.1
Panel 1	027301	2.23	1.89-2.98
Panel 2	124507	12.9	10.1-15.1
Panel 3	125905	34.1	26.3-39.4
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132502	170	129-332
Control 2	132502	3793	2972-6169
Panel 1	027301	31.2	27.7-60.9
Panel 2	124507	117	110-186
Panel 3	125905	2628	2322-4127
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132502	19.7	15.4-28.8
Control 2	132502	440	374-562
Panel 1	027301	10.5	8.38-12.6
Panel 2	124507	38.5	21.8-48.2
Panel 3	125905	320	165-410

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

Review/Approval

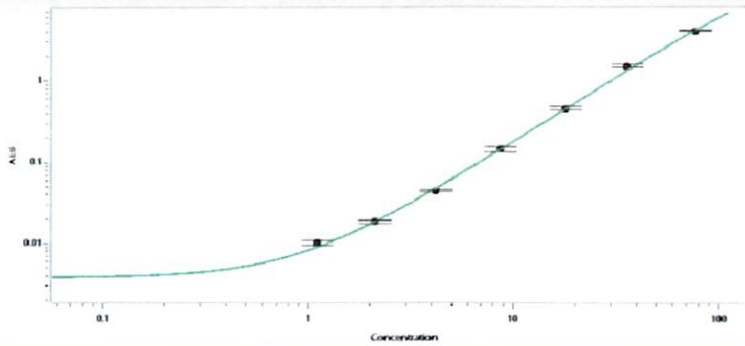
 Director, QC
 29 Sep 2022

Name

Title

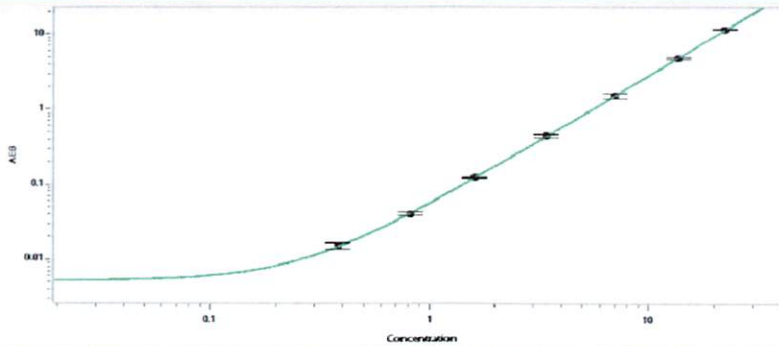
Signature/Date

Abeta 40 Calibration Curve



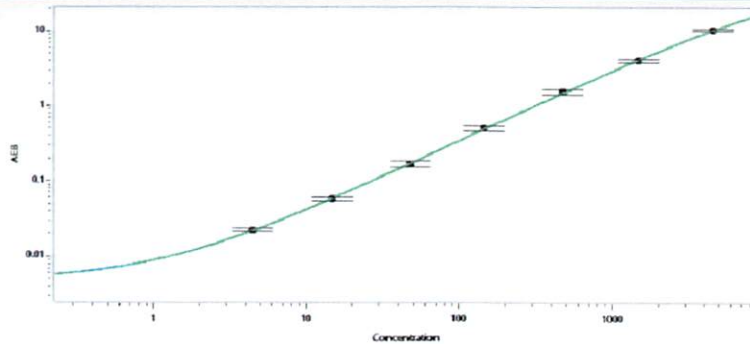
Calibrator Levels (pg/mL)	
A	0.000
B	1.22
C	2.31
D	4.71
E	9.75
F	20.3
G	40.8
H	69.1

Abeta 42 Calibration Curve



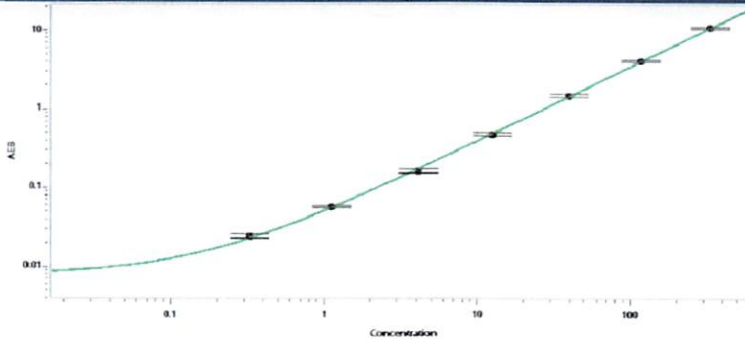
Calibrator Levels (pg/mL)	
A	0.000
B	0.369
C	0.752
D	1.59
E	3.27
F	6.96
G	13.4
H	22.8

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	5.51
C	14.9
D	54.7
E	168
F	599
G	1838
H	4644

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.402
C	1.24
D	4.71
E	14.1
F	49.7
G	147
H	458