

Product Number:	103670
Lot Number:	503859
Expiration:	24-Apr-2024
Platform(s):	HD-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103653	311407	311407A
Detector Reagent	103655	311506	311506A
SBG Reagent	103657	311501	311501A
Plasma Sample Diluent	103659	311504	311504A
CSF Sample Diluent	103727	311505	311505A
Calibrators	103667	315306	315306A
Control 1	103668	315307	315307A
Control 2	103669	315307	315307A
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>



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

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials			
<b>Abeta 40</b>			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315307	12.4	10.2 - 15.3
Control 2	315307	97.4	73.4 - 110
Panel 1	226304	25.1	20.8 - 31.2
Panel 2	225507	56.8	44.5 - 66.8
Panel 3	226908	254	183 - 312
<b>Abeta 42</b>			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315307	4.41	3.60 - 5.40
Control 2	315307	24.9	19.9 - 29.9
Panel 1	226304	2.12	1.65 - 2.97
Panel 2	225507	18.6	16.5 - 24.7
Panel 3	226908	46.6	41.2 - 61.9
<b>GFAP®</b>			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315307	154	122 - 183
Control 2	315307	3083	2415 - 3623
Panel 1	226304	42.6	28.1 - 58.0
Panel 2	225507	264	248 - 372
Panel 3	226908	2172	2078 - 3165
<b>NF-light®</b>			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315307	19.9	15.8 - 23.8
Control 2	315307	478	366 - 549
Panel 1	226304	3.04	1.38 - 4.84
Panel 2	225507	107	85.7 - 128
Panel 3	226908	525	367 - 698

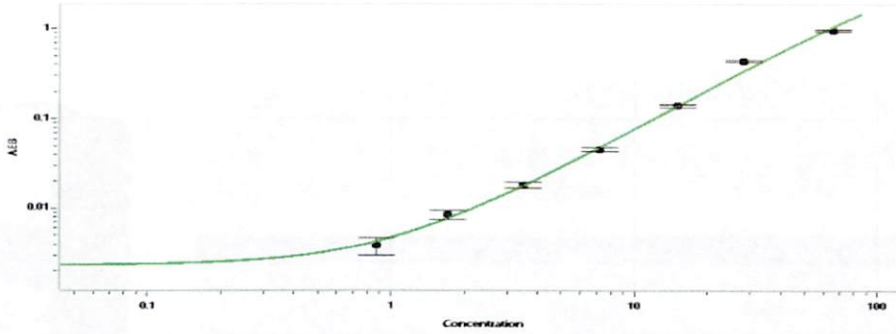
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

ROBERT SOMOGYI QC, MANAGER [Signature] 25 OCT 2023  
 Name Title Signature/Date

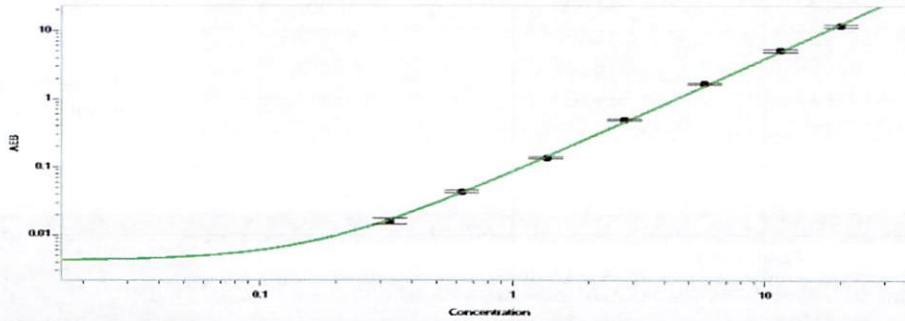


**Abeta 40 Calibration Curve**



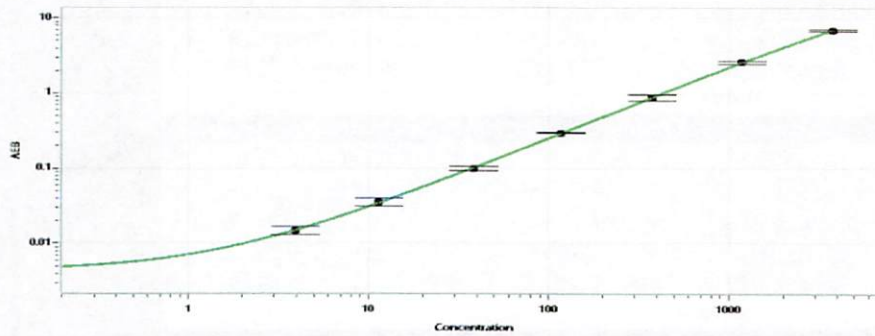
Calibrator Levels (pg/mL)	
A	0.000
B	0.872
C	1.71
D	3.48
E	7.23
F	15.1
G	28.0
H	65.5

**Abeta 42 Calibration Curve**



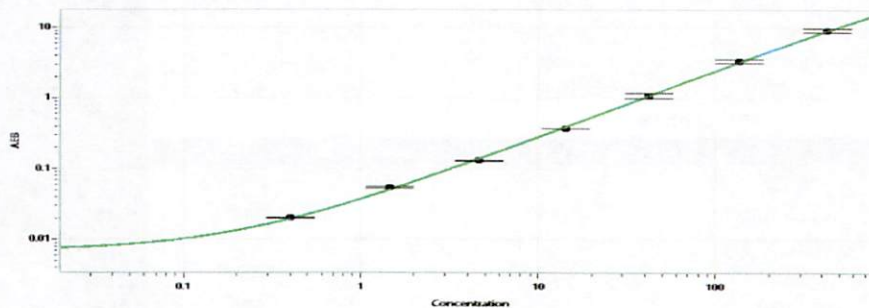
Calibrator Levels (pg/mL)	
A	0.000
B	0.327
C	0.631
D	1.36
E	2.78
F	5.74
G	11.5
H	20.0

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	3.94
C	11.4
D	38.5
E	117
F	375
G	1187
H	3742

**NF-light® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.404
C	1.46
D	4.61
E	14.3
F	41.7
G	135
H	424