

Product Number:	103670
Lot Number:	503504
Expiration:	14-Nov-2023
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	233604
Detector Reagent	103655	233605
SBG Reagent	103657	233609
Plasma Sample Diluent	103659	233610
CSF Sample Diluent	103727	233611
Calibrators	103667	230613
Control 1	103668	230614
Control 2	103669	230614
RGP Reagent	103159	N/A <sup>1</sup>

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

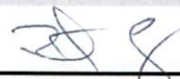
Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	20.1	15.2-22.7
Control 2	230614	133	103-154
Panel 1	226304	26.1	20.3-30.4
Panel 2	225507	58.4	44.3-66.5
Panel 3	226908	245	179-306
Abeta 42			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	5.94	4.65-6.97
Control 2	230614	30.4	24.7-37.0
Panel 1	226304	2.14	1.83-2.83
Panel 2	225507	20.9	16.6-24.9
Panel 3	226908	51.8	41.6-62.4
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	212	164-246
Control 2	230614	4361	3434-5151
Panel 1	226304	43.4	34.0-53.7
Panel 2	225507	319	248-372
Panel 3	226908	2673	2079-3186
NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	19.7	14.9-22.4
Control 2	230614	353	287-430
Panel 1	226304	3.32	2.20-4.28
Panel 2	225507	109	87.1-131
Panel 3	226908	511	402-697

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

Review/Approval

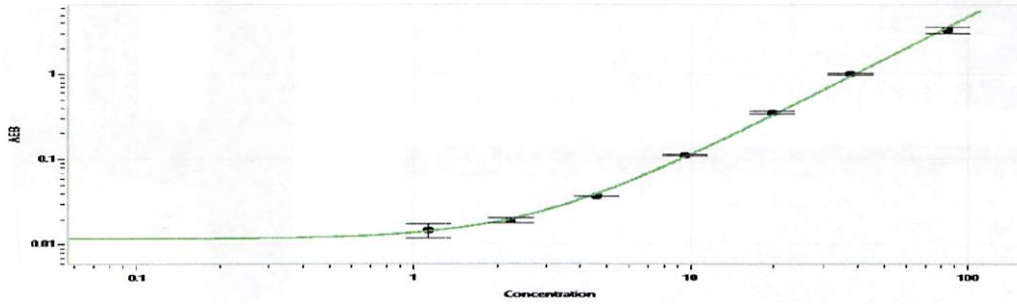
ROBERT SAMOGGI  
Name

QC, MANAGER  
Title

 25052023  
Signature/Date

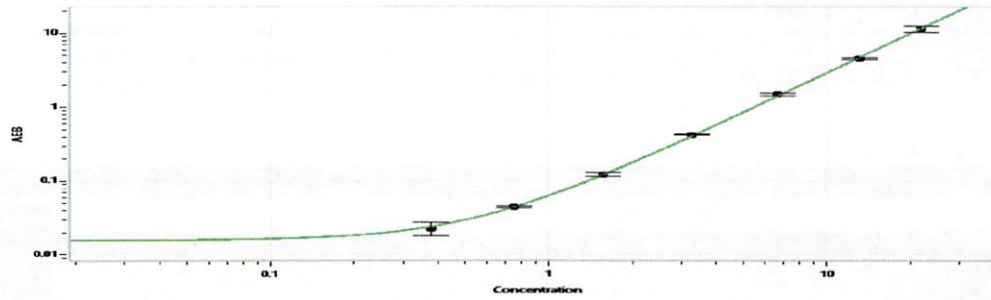


**Abeta 40 Calibration Curve**



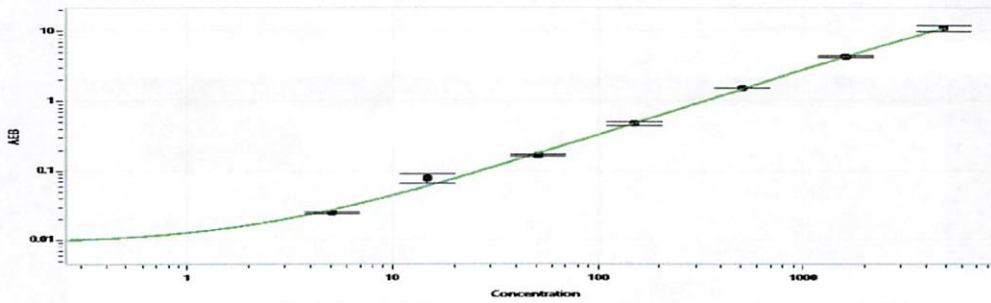
Calibrator Levels (pg/mL)	
A	0.000
B	1.13
C	2.24
D	4.58
E	9.52
F	19.6
G	37.5
H	84.3

**Abeta 42 Calibration Curve**



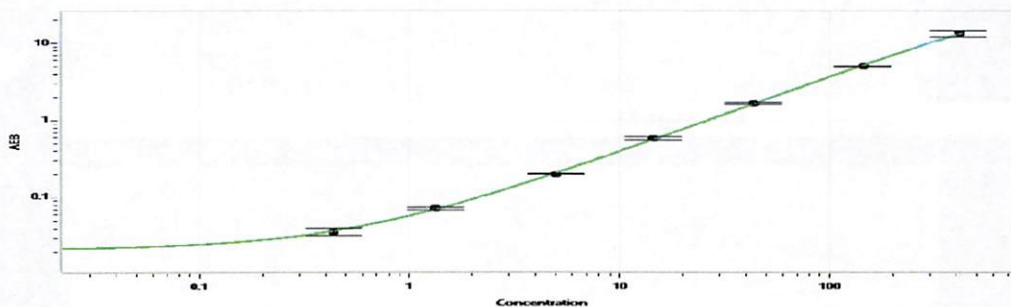
Calibrator Levels (pg/mL)	
A	0.000
B	0.378
C	0.752
D	1.57
E	3.25
F	6.68
G	13.1
H	21.8

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	5.08
C	14.8
D	51.1
E	151
F	508
G	1615
H	4821

**NF-light® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.438
C	1.33
D	4.96
E	14.4
F	43.2
G	144
H	411