

Product Number:	103670
Lot Number:	503420
Expiration:	9-Jun-2023
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	230512
Detector Reagent	103655	230612
SBG Reagent	103657	221004
Plasma Sample Diluent	103659	216004
CSF Sample Diluent	103727	216005
Calibrators	103667	230613
Control 1	103668	230614
Control 2	103669	230614
RGP Reagent	103159	N/A ¹

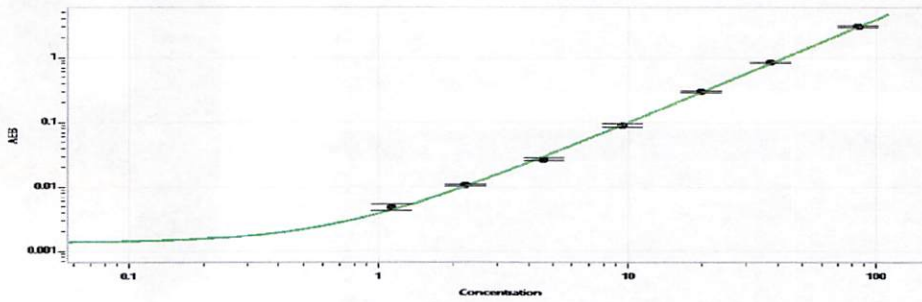
¹ RGP is not Kit Lot Specific

Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	18.3	14.8-22.2
Control 2	230614	134	103-155
Panel 1	027301	25.3	20.8-36.1
Panel 2	225507	58.4	43.7-65.6
Panel 3	226908	264	172-313
Abeta 42			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	5.73	4.60-6.90
Control 2	230614	31.3	24.9-37.3
Panel 1	027301	2.17	1.89-2.99
Panel 2	225507	20.4	16.6-24.9
Panel 3	226908	52.3	41.6-62.5
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	210	164-245
Control 2	230614	4914	3541-5320
Panel 1	027301	47.8	29.7-59.6
Panel 2	225507	331	246-369
Panel 3	226908	2837	2049-3197
NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	18.8	14.7-22.1
Control 2	230614	392	295-442
Panel 1	027301	10.5	8.38-12.6
Panel 2	225507	109	87.1-131
Panel 3	226908	550	407-711

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

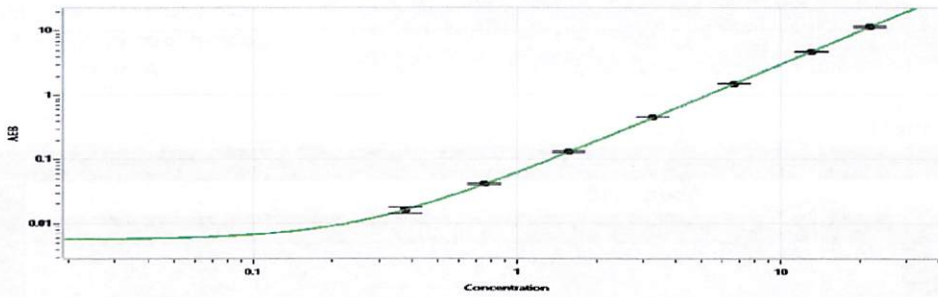
Review/Approval: Sonia Khillan Name QC Manager Title Sonia Khillan | 28 Nov 2022 Signature/Date

Abeta 40 Calibration Curve



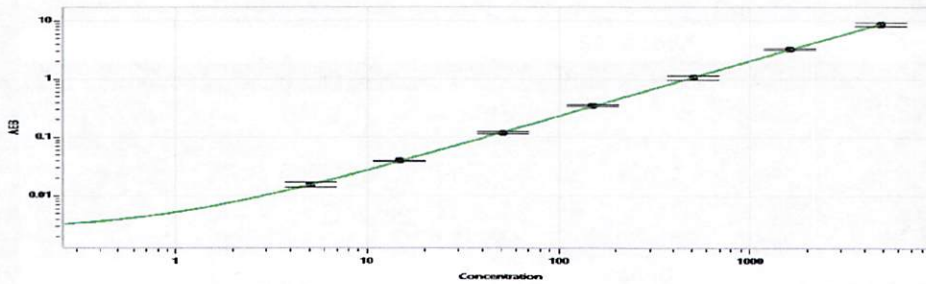
Calibrator Levels (pg/mL)	
A	0.000
B	1.13
C	2.24
D	4.58
E	9.52
F	19.6
G	37.5
H	84.3

Abeta 42 Calibration Curve



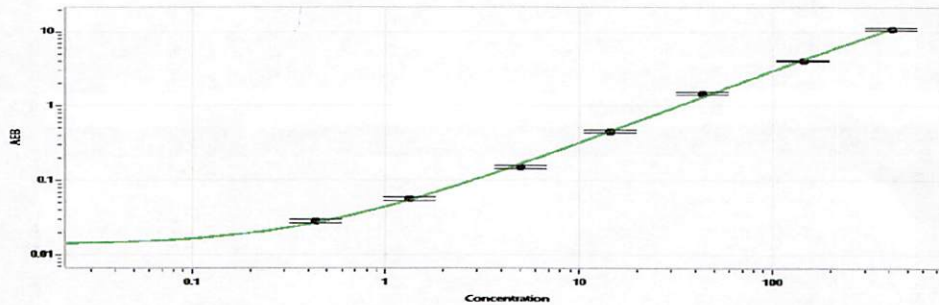
Calibrator Levels (pg/mL)	
A	0.000
B	0.378
C	0.752
D	1.57
E	3.25
F	6.68
G	13.1
H	21.8

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