

Product Number:	103670
Lot Number:	503745
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	232510
Detector Reagent	103655	232703
SBG Reagent	103657	232704
Plasma Sample Diluent	103659	232705
CSF Sample Diluent	103727	233306
Calibrators	103667	230613
Control 1	103668	230614
Control 2	103669	230614
RGP Reagent	103159	N/A ¹

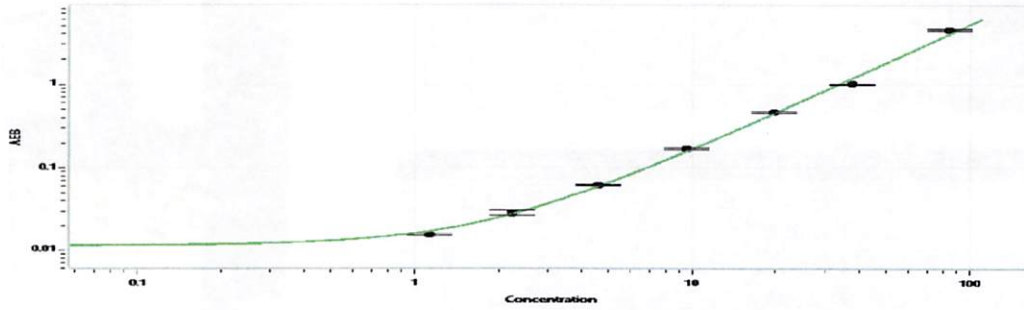
¹ RGP is not Kit Lot Specific

Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	17.7	15.0-22.4
Control 2	230614	119	101-152
Panel 1	226304	26.5	20.4-30.7
Panel 2	225507	57.1	44.5-66.8
Panel 3	226908	273	183-312
Abeta 42			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	5.30	4.56-6.84
Control 2	230614	28.5	24.3-36.5
Panel 1	226304	1.96	1.65-2.86
Panel 2	225507	19.6	16.5-24.7
Panel 3	226908	49.3	41.2-61.9
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	215	166-248
Control 2	230614	4257	3428-5142
Panel 1	226304	35.7	30.1-54.3
Panel 2	225507	309	248-372
Panel 3	226908	2563	2078-3165
NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	230614	17.0	14.6-22.0
Control 2	230614	318	280-421
Panel 1	226304	1.96	1.35-4.62
Panel 2	225507	98.2	85.7-128
Panel 3	226908	450	367-698

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

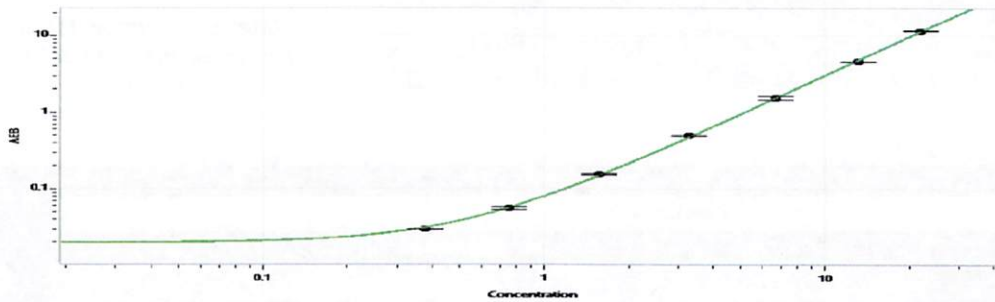
Review/Approval	<u>ROBERT SONOGTI</u>	<u>QC, MANAGER</u>	<u>[Signature]</u> 25 Oct 2023
	Name	Title	Signature/Date

Abeta 40 Calibration Curve



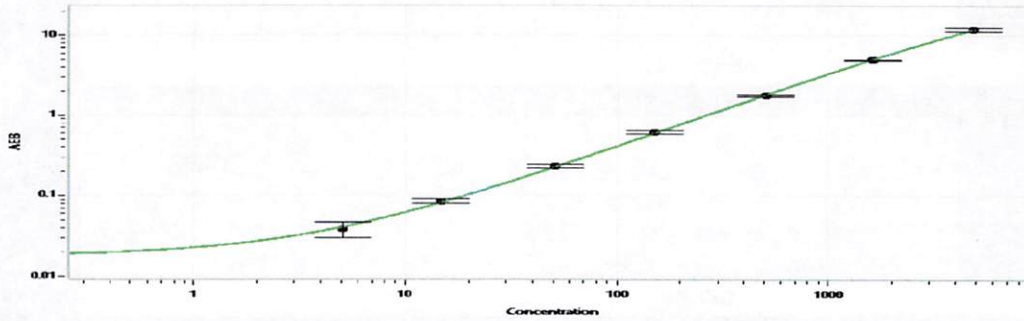
Calibrator Levels (pg/mL)	
A	0.000
B	1.13
C	2.24
D	4.58
E	9.52
F	19.6
G	37.5
H	84.3

Abeta 42 Calibration Curve



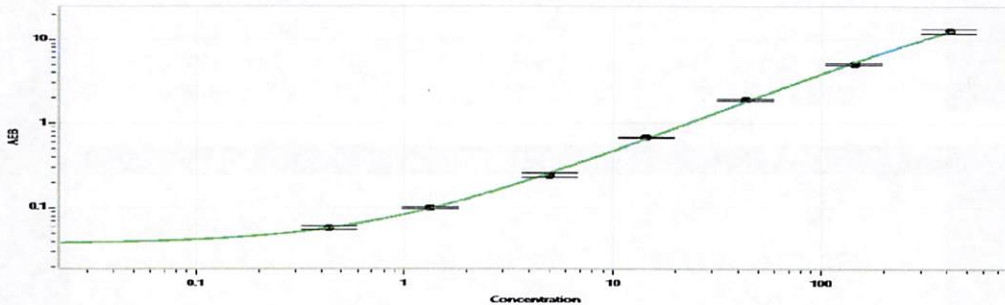
Calibrator Levels (pg/mL)	
A	0.000
B	0.378
C	0.752
D	1.57
E	3.25
F	6.68
G	13.1
H	21.8

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