

Product Number:	103670
Lot Number:	503768
Expiration:	11-Jan-2024
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	301108
Detector Reagent	103655	301208
SBG Reagent	103657	301717
Plasma Sample Diluent	103659	301304
CSF Sample Diluent	103727	301713
Calibrators	103667	304107
Control 1	103668	304108
Control 2	103669	304108
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	304108	17.7	14.2-21.3
Control 2	304108	120	96.2-144
Panel 1	226304	28.1	20.8-31.2
Panel 2	225507	59.3	44.5-66.8
Panel 3	226908	275	183-312
Abeta 42			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	304108	5.86	4.55-6.82
Control 2	304108	30.4	24.1-36.2
Panel 1	226304	2.56	1.65-2.97
Panel 2	225507	22.2	16.5-24.7
Panel 3	226908	52.6	41.2-61.9
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	304108	185	147-220
Control 2	304108	3643	2814-4221
Panel 1	226304	47.2	28.1 - 58.0
Panel 2	225507	311	248-372
Panel 3	226908	2340	2078-3165
NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	304108	23.7	18.5-27.7
Control 2	304108	483	345-543
Panel 1	226304	3.73	1.38-4.84
Panel 2	225507	115	85.7-128
Panel 3	226908	493	367-698

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

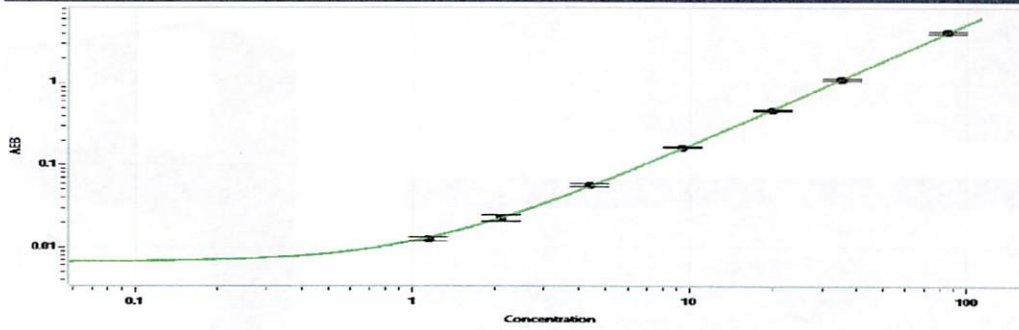
Review/Approval

ROBERT SONOGTI
Name

QC, MANAGER
Title

[Signature] 25 Oct 2023
Signature/Date

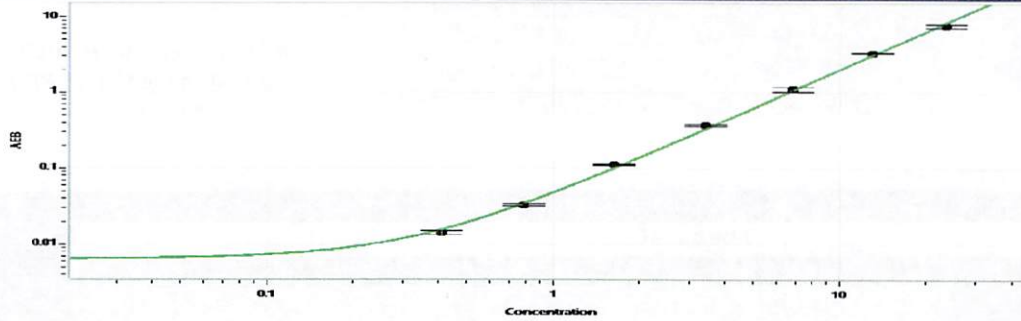
Abeta 40 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.15
C	2.10
D	4.38
E	9.40
F	19.8
G	35.2
H	85.2

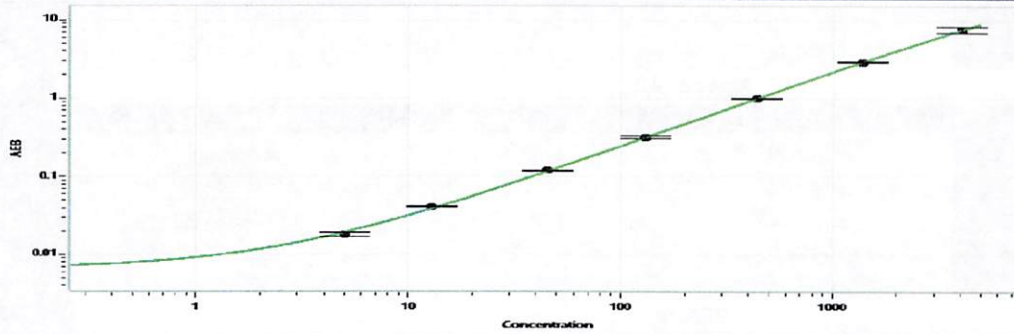
Abeta 42 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.407
C	0.796
D	1.63
E	3.41
F	6.89
G	13.1
H	23.6

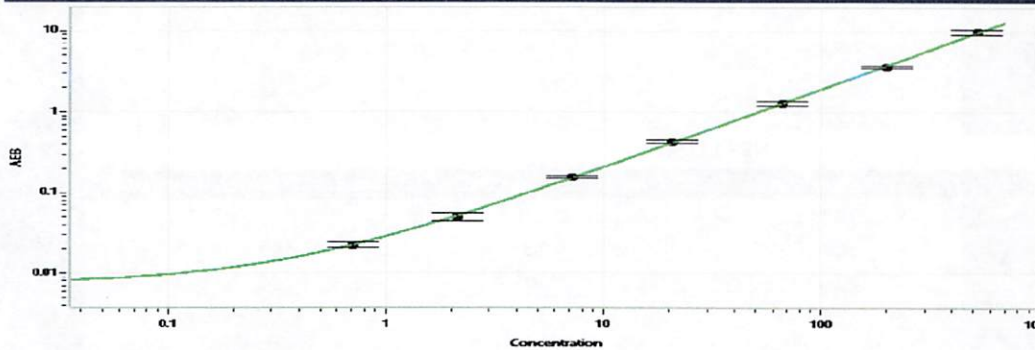
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	5.02
C	12.8
D	45.3
E	132
F	441
G	1390
H	4064

NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.706
C	2.13
D	7.17
E	20.7
F	66.4
G	200
H	518