

Product Number:	103670
Lot Number:	503246
Expiration:	6-Nov-2022
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	131001
Detector Reagent	103655	133405
SBG Reagent	103657	133411
Plasma Sample Diluent	103659	133515
CSF Sample Diluent	103727	133516
Calibrators	103667	134101
Control 1	103668	134102
Control 2	103669	134102
RGP Reagent	103159	N/A

¹ RGP is not Kit Lot Specific

Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134102	18.1	7.10-27.3
Control 2	134102	124	91.0-136
Panel 1	027301	25.0	17.4-36.4
Panel 2	124507	84.9	69.2-104
Panel 3	125905	116	92.1-138
Abeta 42			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134102	5.97	4.76-7.13
Control 2	134102	32.6	26.2-39.4
Panel 1	027301	1.90	1.87-3.00
Panel 2	124507	11.8	10.1-15.1
Panel 3	125905	30.6	26.3-39.4
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134102	221	168-251
Control 2	134102	4082	3270-4905
Panel 1	027301	44.6	32.9-57.3
Panel 2	124507	156	110-186
Panel 3	125905	2856	2322-4127
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134102	17.8	13.4-20.1
Control 2	134102	341	304-455
Panel 1	027301	9.67	7.13-12.9
Panel 2	124507	33.6	21.8-48.2
Panel 3	125905	251	165-410

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

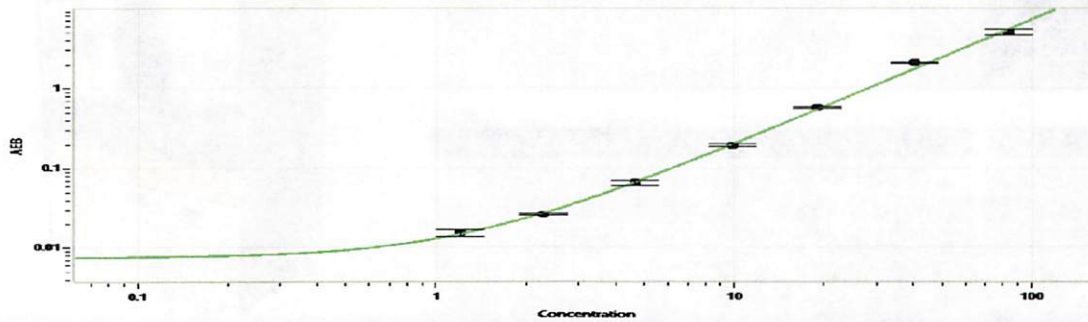
Review/Approval

Wenkui Ts. Name

QC Supervisor Title

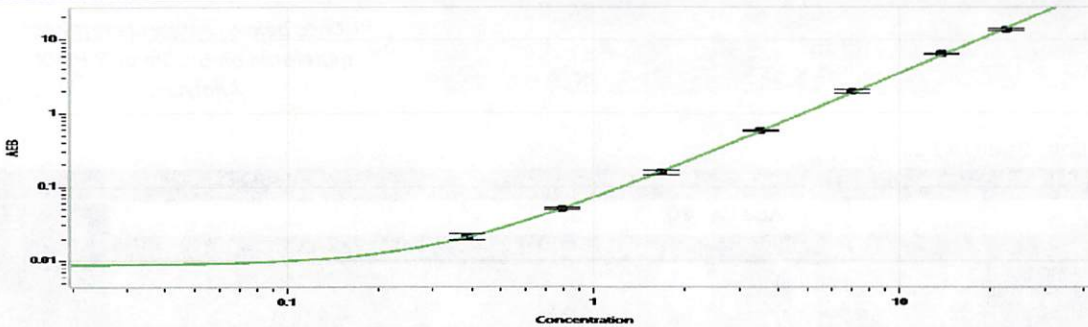
[Signature] 14 Jun 2022 Signature/Date

Abeta 40 Calibration Curve



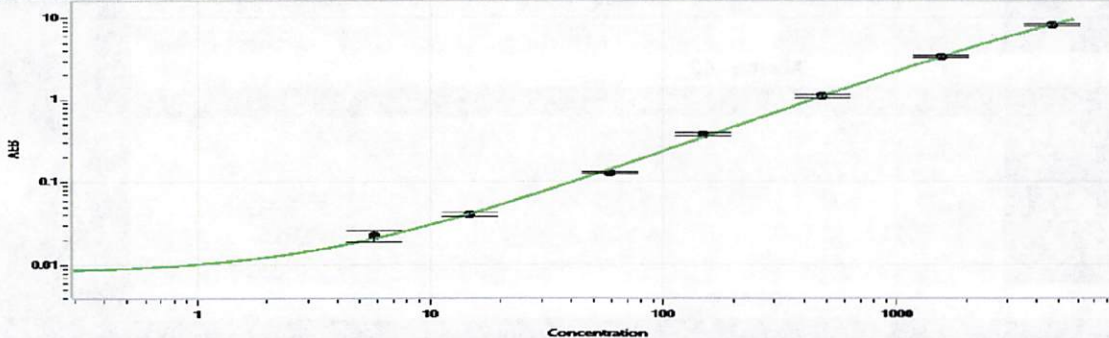
Calibrator Levels (pg/mL)	
A	0.000
B	1.21
C	2.31
D	4.62
E	9.81
F	19.1
G	40.4
H	84.0

Abeta 42 Calibration Curve



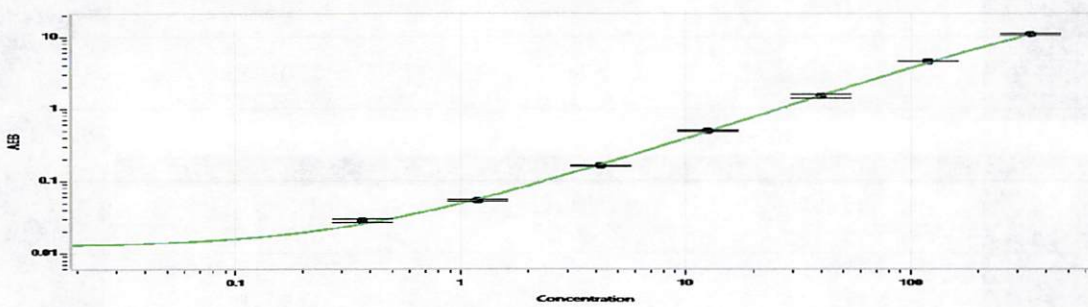
Calibrator Levels (pg/mL)	
A	0.000
B	0.388
C	0.787
D	1.66
E	3.51
F	7.04
G	13.8
H	22.4

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	5.77
C	14.8
D	58.6
E	149
F	477
G	1552
H	4646

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.370
C	1.19
D	4.18
E	12.5
F	40.1
G	120
H	344