

Product Number:	103670
Lot Number:	503236
Expiration:	18-Sep-2022
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	126101
Detector Reagent	103655	128801
SBG Reagent	103657	126105
Plasma Sample Diluent	103659	126202
CSF Sample Diluent	103727	126203
Calibrators	103667	128603
Control 1	103668	128604
Control 2	103669	128604
RGP Reagent	103159	N/A

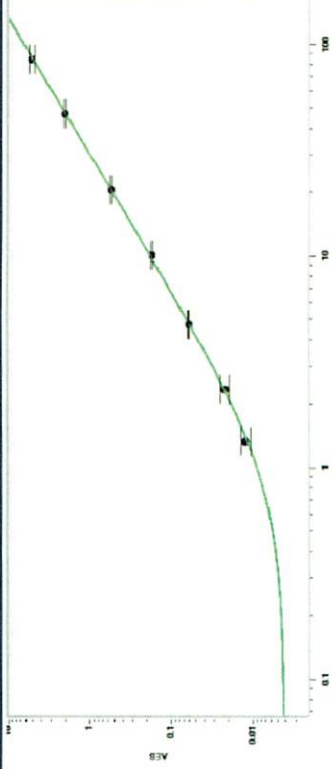
¹ RGP is not Kit Lot Specific

Release Materials			
Abeta 40			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	128604	17.8	14.2-21.3
Control 2	128604	130	106-159
Panel 1	027301	23.6	17.4-36.4
Panel 2	124507	89.3	69.2-104
Panel 3	125905	118	92.1-138
Abeta 42			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	128604	5.87	4.59-6.88
Control 2	128604	31.3	24.9-37.3
Panel 1	027301	2.08	1.87-3.00
Panel 2	124507	12.3	10.1-15.1
Panel 3	125905	31.4	26.3-39.4
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	128604	234	177-266
Control 2	128604	4154	3309-4963
Panel 1	027301	49.0	32.9-57.3
Panel 2	124507	165	110-186
Panel 3	125905	3137	2322-4127
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	128604	21.2	16.6-24.9
Control 2	128604	395	317-476
Panel 1	027301	9.28	7.13-12.9
Panel 2	124507	31.6	21.8-48.2
Panel 3	125905	232	165-410

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

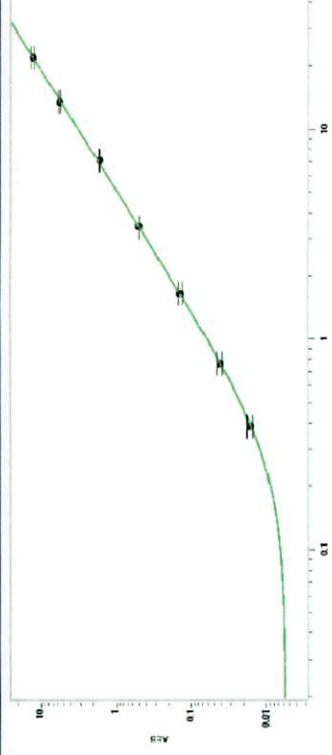
Review/Approval	<u>Shivani Goel</u>	<u>QC</u>	<u>Shivani Goel 17 Mar 2022</u>
	Name	Title	Signature/Date

Abeta 40 Calibration Curve



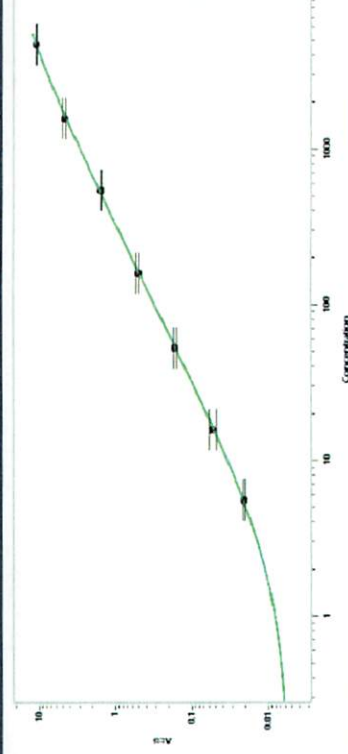
Calibrator Levels (pg/ml)
A 0.000
B 1.33
C 2.32
D 4.73
E 10.1
F 20.5
G 46.5
H 83.8

Abeta 42 Calibration Curve



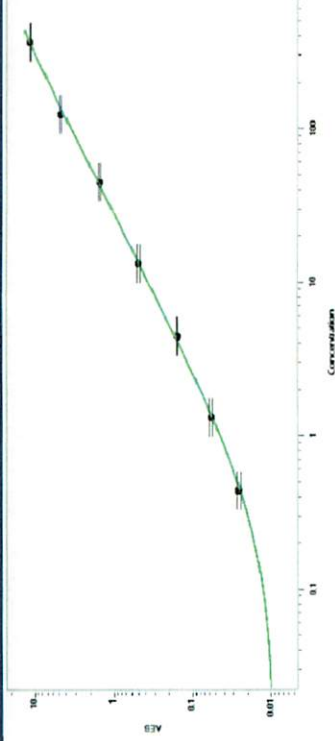
Calibrator Levels (pg/ml)
A 0.000
B 0.383
C 0.764
D 1.64
E 3.41
F 7.03
G 13.3
H 21.7

GFAP® Calibration Curve



Calibrator Levels (pg/ml)
A 0.000
B 5.54
C 15.7
D 52.8
E 158
F 543
G 1559
H 4683

NF-light® Calibration Curve



Calibrator Levels (pg/ml)
A 0.000
B 0.441
C 1.31
D 4.39
E 13.3
F 44.3
G 124
H 366