

Product Number:	104465
Lot Number:	504182
Expiration:	23-Aug-2025
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104448	405409
Detector Reagent	104450	419710
SBG Reagent	104452	407306
Calibrator Diluent Reagent	104453	405911
Plasma Sample Diluent Reagent	104647	406008
CSF Sample Diluent Reagent	104648	406009
Control Diluent Reagent	104649	405912
Calibrator Concentrate	104466	402917
Control 1 Concentrate	104468	402918
Control 2 Concentrate	104469	402919
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 07 Aug 2024

Signature/Date

Release Materials*

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402918	4.10	2.76 - 5.13
Control 2	402919	60.6	39.3 - 72.9

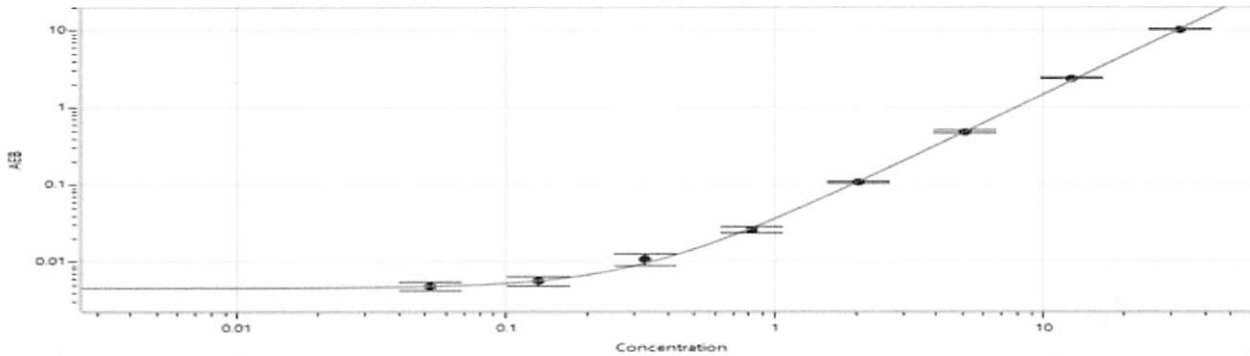
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibrator Levels (pg/mL)

A	0.000
B	0.052
C	0.131
D	0.329
E	0.821
F	2.05
G	5.13
H	12.8
I	32.1

Calibration Curve



Release Materials*

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402918	2.30	1.42 - 2.64
Control 2	402919	32.2	21.1 - 39.2

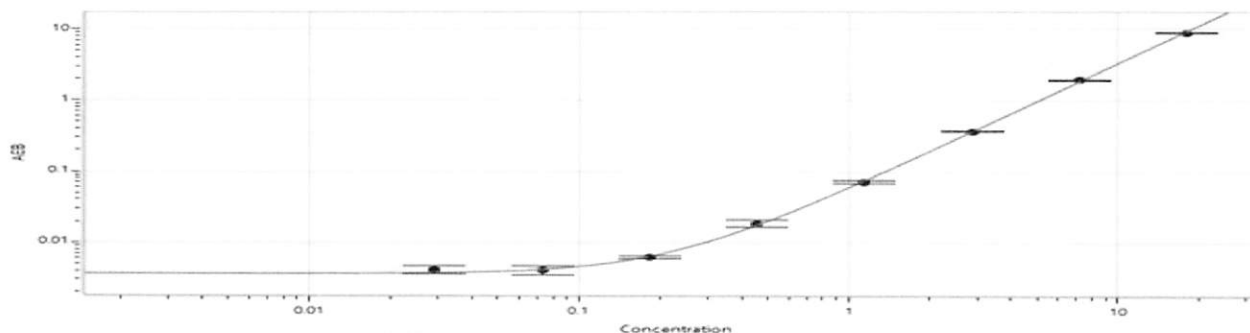
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibrator Levels (pg/mL)

A	0.000
B	0.029
C	0.073
D	0.183
E	0.457
F	1.14
G	2.86
H	7.14
I	17.8

Calibration Curve



Release Materials*

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402918	95.8	65.9 - 122
Control 2	402919	1320	832 - 1545

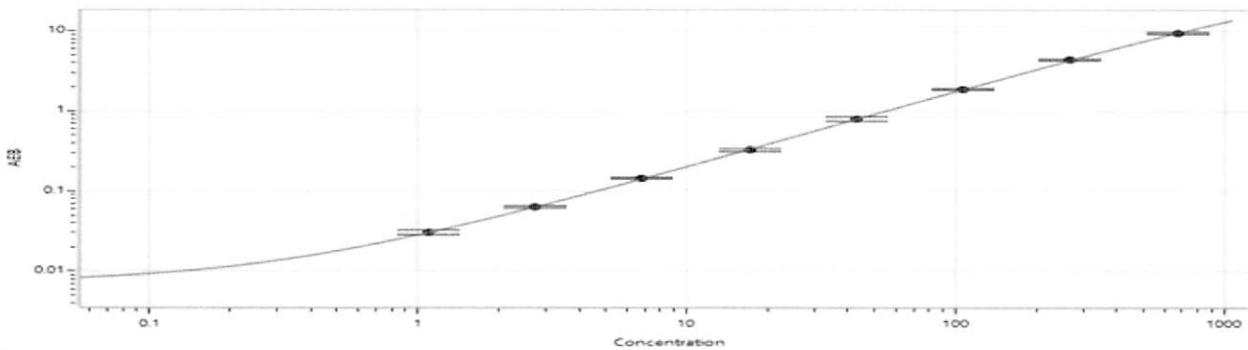
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibrator Levels (pg/mL)

A	0.000
B	1.10
C	2.75
D	6.86
E	17.2
F	42.8
G	107
H	268
I	670

Calibration Curve



Release Materials*

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402918	92.7	54.5 - 101
Control 2	402919	1314	713 - 1324

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibrator Levels (pg/mL)

A	0.000
B	0.866
C	2.17
D	5.41
E	13.5
F	33.8
G	84.6
H	211
I	529

Calibration Curve

