

Product Number:	103670
Lot Number:	502542
Expiration:	14-Jul-2021
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103653	026603
Detector Reagent	103655	026807
SBG Reagent	103657	026711
Plasma Sample Diluent	103659	026708
CSF Sample Diluent	103727	026802
Calibrators	103667	028102
Control 1	103668	028103
Control 2	103669	028103
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0154).

**Release Materials**

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028103	15.5	12.6-18.9
Control 2	028103	118	92.5-139
Panel 1	027301	26.5	23.3-35.0
Panel 2	018109	53.0	43.6-65.4
Panel 3	018109	137	109-164

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028103	5.82	4.66-7.00
Control 2	028103	32.3	25.6-38.3
Panel 1	027301	2.30	2.01-3.03
Panel 2	018109	20.2	16.8-25.2
Panel 3	018109	45.9	36.2-54.3

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028103	235	183-275
Control 2	028103	5055	3761-5641
Panel 1	027301	44.9	37.8-56.7
Panel 2	018109	547	446-669
Panel 3	018109	2459	1892-2940

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028103	33.1	20.6-37.0
Control 2	028103	669	393-774
Panel 1	027301	10.82	8.37-12.6
Panel 2	018109	77.4	59.7-89.5
Panel 3	018109	415	305-473

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

Review/Approval

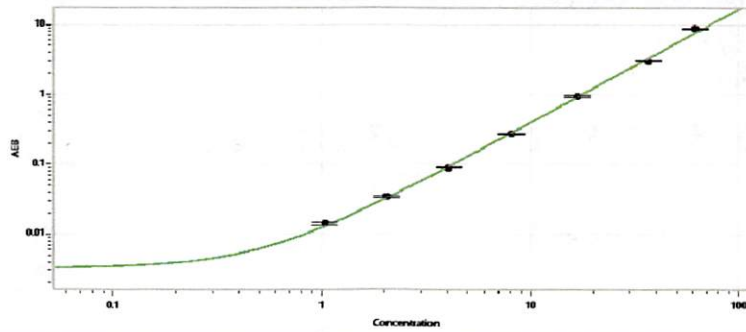
Shivani Goel  
Name

QC Manager  
Title

Shivani Goel Boct2020  
Signature/Date

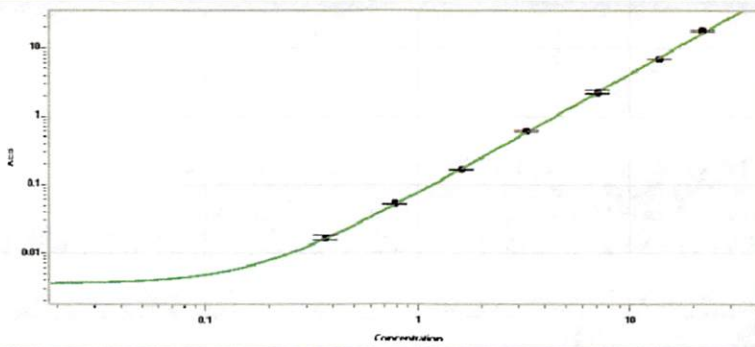


**Abeta 40 Calibration Curve**



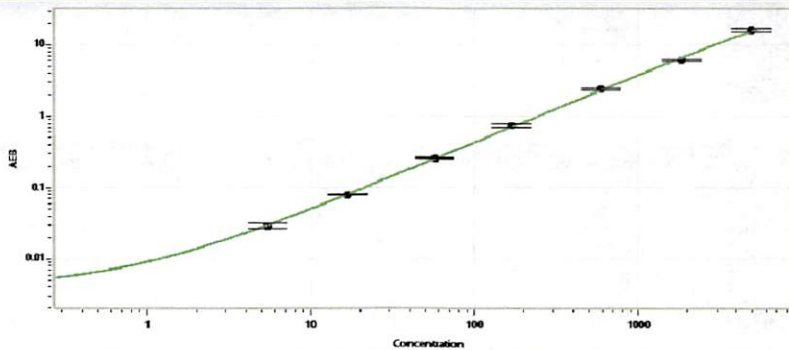
Calibrator Levels (pg/mL)	
A	0.000
B	1.04
C	2.05
D	4.09
E	8.14
F	16.8
G	36.6
H	61.4

**Abeta 42 Calibration Curve**



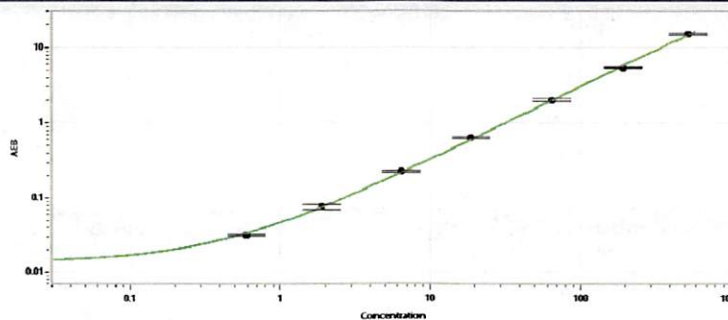
Calibrator Levels (pg/mL)	
A	0.000
B	0.367
C	0.782
D	1.62
E	3.25
F	7.07
G	13.8
H	22.0

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	5.41
C	16.7
D	56.9
E	166
F	589
G	1827
H	4920

**NF-light® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.597
C	1.88
D	6.42
E	18.7
F	64.4
G	195
H	528