

Product Number:	104797
Lot Number:	504275
Expiration:	23-Nov-2025
Platform(s):	HD-X



Component	Part Number	Lot Number
Bead Reagent	104778	414404
Detector Reagent	104780	414910
SBG Reagent	104782	414911
Plasma Sample Diluent	104783	414512
CSF Sample Diluent	104784	414411
Calibrator Diluent	104786	414412
Control Diluent	104787	414413
Calibrator Concentrate	104793	409512
Control 1 Concentrate	104795	409513
Control 2 Concentrate	104796	409514
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

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

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	409513	7.12	5.62 - 10.4
Control 2 Concentrate	409514	318	225 - 418
Panel 1	410108	2.32	1.80 - 3.35
Panel 2	410108	21.7	16.6 - 30.8
Panel 3	410108	139	111 - 206
Panel 4	410108	2.04	1.48 - 2.75
Panel 5	410108	24.3	17.9 - 33.2
Panel 6	410108	159	120 - 223

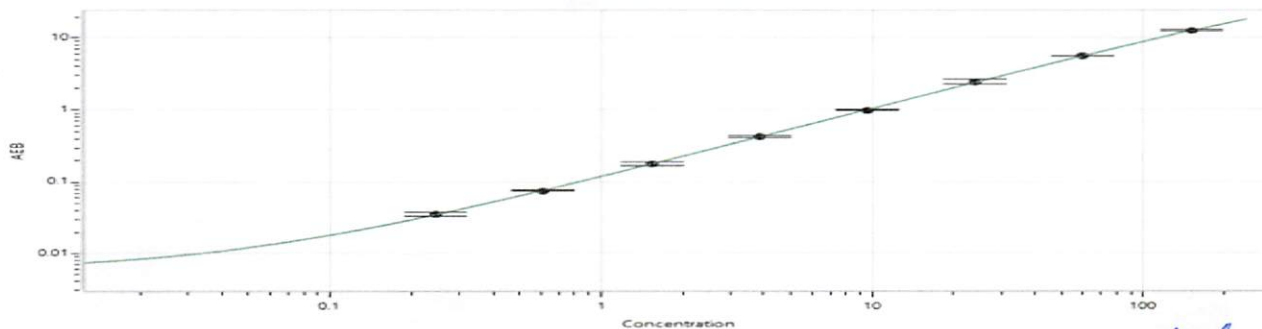
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1465
Control 1 Concentrate Reconstitution Volume (uL)	1465
Control 2 Concentrate Reconstitution Volume (uL)	1465

Calibrator Levels (pg/mL)

A	0
B	0.246
C	0.614
D	1.54
E	3.84
F	9.60
G	24.0
H	60.0
I	150

Calibration Curve



Review/Approval

Cody Carlson
Name Title

[Signature] 18 JUN 2024
Signature/Date