

Product Number:	104691
Lot Number:	504476
Expiration:	13-Feb-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104752	501712	501712A
Detector Reagent	104675	502703	502703A
SBG Reagent	104677	502704	502704A
Plasma Sample Diluent	104719	502114	502114A
CSF Sample Diluent	104720	502115	502115A
Calibrator Diluent	104678	422628	N/A
Control Diluent	104718	422629	N/A
Calibrator Concentrate	104687	402944	N/A
Control 1 Concentrate	104689	402945	N/A
Control 2 Concentrate	104690	402946	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

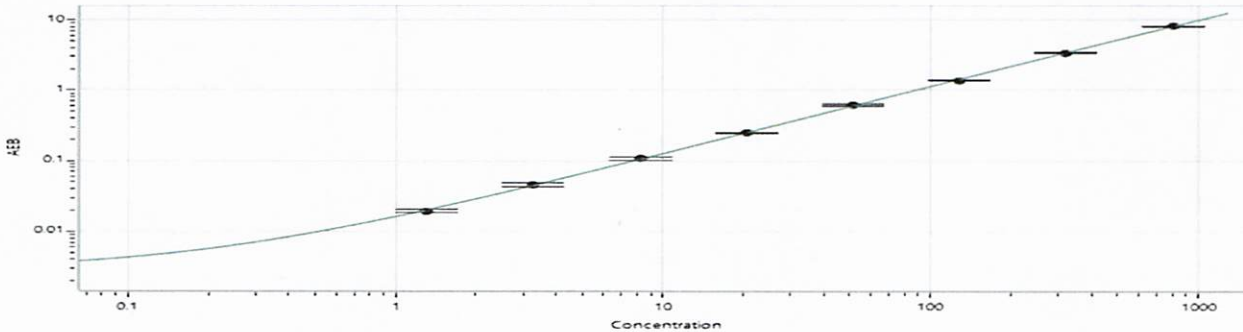
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	402945	96.9	78.6 - 146
Control 2 Concentrate	402946	1262	964 - 1790

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

Calibrator Concentrate Reconstitution Volume (uL)	1053
Control 1 Concentrate Reconstitution Volume (uL)	1053
Control 2 Concentrate Reconstitution Volume (uL)	1053

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.31
C	3.28
D	8.19
E	20.5
F	51.2
G	128
H	320
I	800

Review/Approval

Kyle Giroux
Name

Senior Document Control Specialist, QA
Title

Kyle Giroux 05 Mar 2025
Signature/Date