

Product Number:	104691
Lot Number:	504630
Expiration:	23-Nov-2026
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104752	516802
Detector Reagent	104675	524708
SBG Reagent	104677	524709
Plasma Sample Diluent	104719	514709
CSF Sample Diluent	104720	514315
Calibrator Diluent	104678	516805
Control Diluent	104718	516806
Calibrator Concentrate	104687	403918
Control 1 Concentrate	104689	403919
Control 2 Concentrate	104690	403920
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



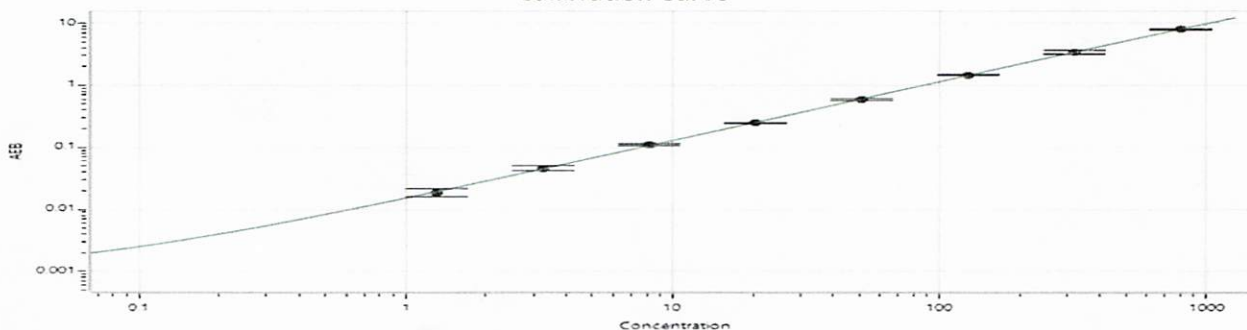
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	403919	99.9	74.6 - 139
Control 2 Concentrate	403920	1497	987 - 1833

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

Calibrator Concentrate Reconstitution Volume (uL)	1393
Control 1 Concentrate Reconstitution Volume (uL)	1393
Control 2 Concentrate Reconstitution Volume (uL)	1393

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

Dipesh Patel Sr. QA Specialist
Name Title

AT 30 Sep 2025
Signature/Date