

Product Number:	104570
Lot Number:	504654
Expiration:	2-Jan-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	518407
Detector Reagent	104548	519001
SBG Reagent	104550	521905
Sample Diluent	104552	518308
Calibrator Diluent Reagent	104745	518309
Control Diluent Reagent	104554	518310
Calibrator Concentrate	104566	406408
Control 1 Concentrate	104568	406409
Control 2 Concentrate	104569	406410
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.

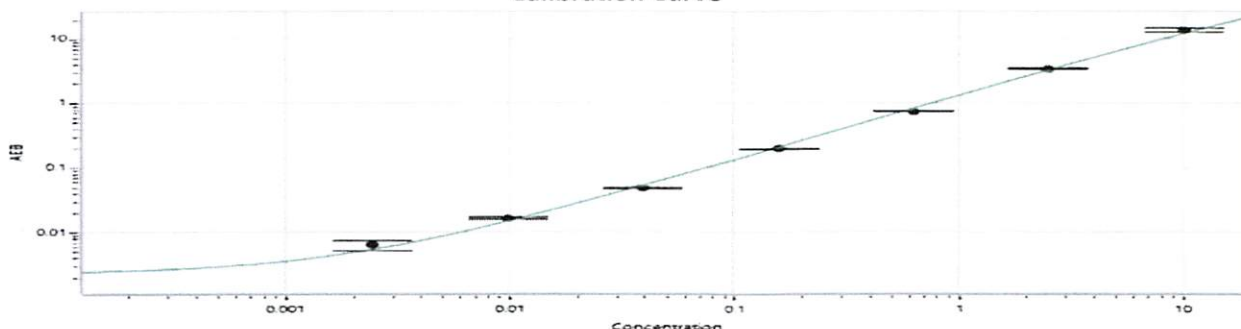
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	406409	0.758	0.517 - 0.961
Control 2 Concentrate	406410	8.06	5.15 - 9.56

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

Calibrator Concentrate Reconstitution Volume (uL)	729
Control 1 Concentrate Reconstitution Volume (uL)	729
Control 2 Concentrate Reconstitution Volume (uL)	729

Calibrator Levels (pg/mL)	
---------------------------	--

Calibration Curve



A	0.000
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0

Review/Approval

Dhruvi Patel QA Specialist II, QA *DP* 11 Sep 2025
 Name Title Signature/Date