

Product Number:	104570
Lot Number:	504657
Expiration:	2-Jan-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	521810
Detector Reagent	104548	523309
SBG Reagent	104550	523310
Sample Diluent	104552	518315
Calibrator Diluent Reagent	104745	518317
Control Diluent Reagent	104554	518316
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

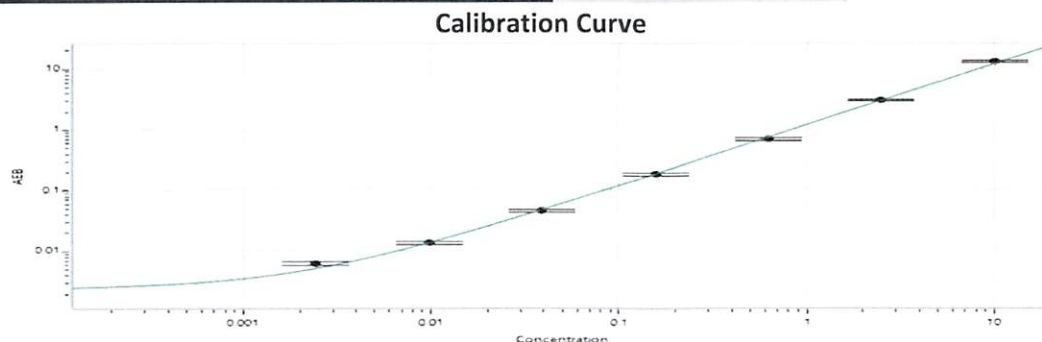
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	405211	0.787	0.494 - 0.917
Control 2 Concentrate	405212	8.13	4.91 - 9.12

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

Calibrator Concentrate Reconstitution Volume (uL)	826
Control 1 Concentrate Reconstitution Volume (uL)	826
Control 2 Concentrate Reconstitution Volume (uL)	826

**Calibrator Levels
(pg/mL)**



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

Dhwani Patel QA Specialist II, QA DP | 19 Sep 2025
Name Title Signature/Date