

Product Number:	104364
Lot Number:	504231
Expiration:	7-Aug-2026
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	104352	416217
Detector Reagent	104353	416212
SBG Reagent	104354	420509
Sample Diluent Reagent	104355	416201
Calibrator Diluent Reagent	104358	415913
Control Diluent Reagent	104359	415914
Calibrator Concentrate	104361	404519
Control 1 Concentrate	104362	404520
Control 2 Concentrate	104363	404521
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.

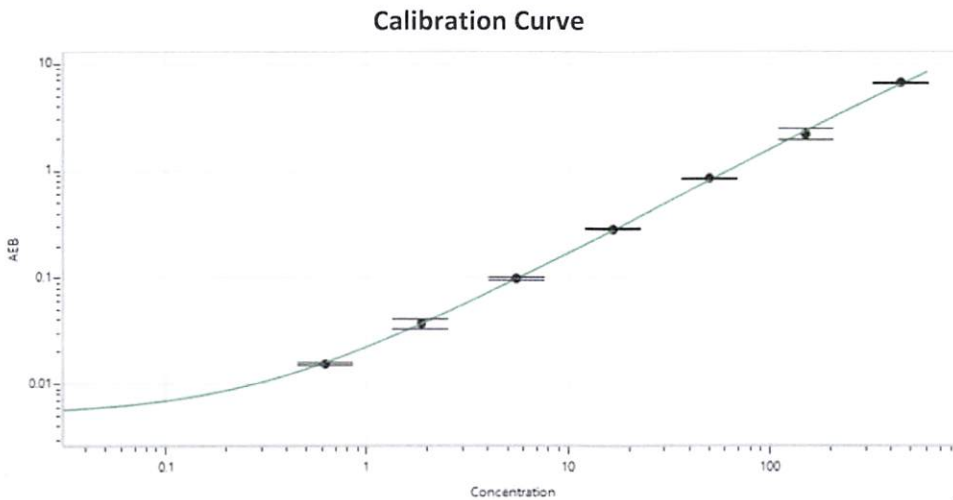
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	404520	4.24	2.67 - 7.20
Control 2	404521	343	250 - 588

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

Calibrator Concentrate Reconstitution Volume (uL)	1480
Control 1 Concentrate Reconstitution Volume (uL)	1480
Control 2 Concentrate Reconstitution Volume (uL)	1480

Calibrator Levels (pg/mL)	
A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450



Signature: [Handwritten Signature] Date: 19 AUG 2024

Review/Approval: Robert Somogyi, QC Manager