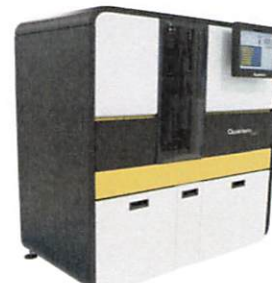


Product Number:	104364
Lot Number:	504465
Expiration:	15-Jun-2026
Platform(s):	HD-X, SR-X

Component	Part Number	Lot Number
Bead Reagent	104352	433001
Detector Reagent	104353	410702
SBG Reagent	104354	433806
Sample Diluent Reagent	104355	410704
Calibrator Diluent Reagent	104358	410614
Control Diluent Reagent	104359	410615
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.

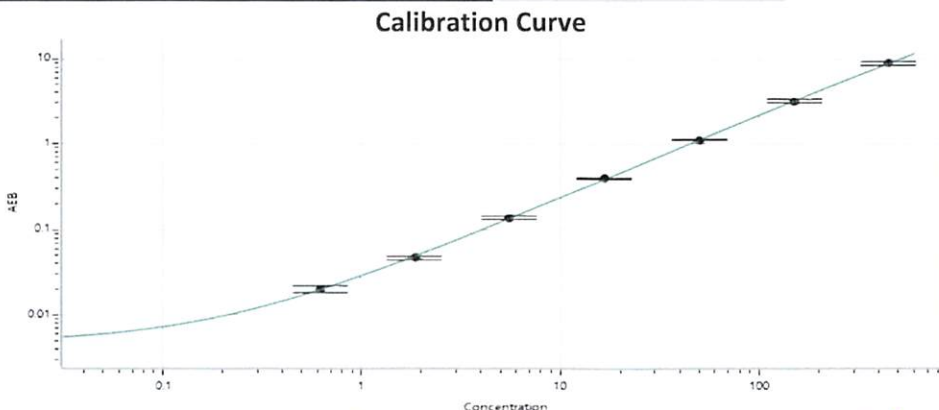
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	404520	3.74	1.68 - 6.64
Control 2	404521	338	275 - 511

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

Calibrator Concentrate Reconstitution Volume (uL)	1323
Control 1 Concentrate Reconstitution Volume (uL)	1323
Control 2 Concentrate Reconstitution Volume (uL)	1323

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



Review/Approval

TAM LEE DOC CONTROL II, QA *Tam Lee* 12 DEC 2024

Name

Title

Signature/Date