

Product Number:	104955
Lot Number:	504427
Expiration:	16-Mar-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104935	428305
Detector Reagent	104937	431301
SBG Reagent	104939	431302
Sample Diluent	104941	426815
Calibrator Diluent	104944	426014
Control Diluent	104945	426015
Calibrator Concentrate	104951	416213
Control 1 Concentrate	104953	416215
Control 2 Concentrate	104954	416216
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

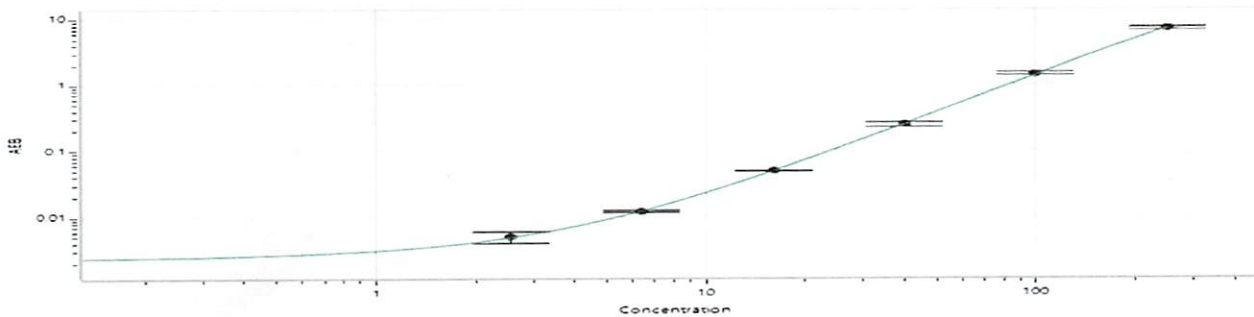
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	416215	44.1	35.8 - 66.5
Control 2 Concentrate	416216	394	276 - 513

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

Calibrator Concentrate Reconstitution Volume (uL)	940
Control 1 Concentrate Reconstitution Volume (uL)	940
Control 2 Concentrate Reconstitution Volume (uL)	940

Calibration Curve



Calibrator Levels (pg/mL)

A	0
B	2.56
C	6.40
D	16.0
E	40.0
F	100
G	250

Review/Approval

Dipesh Patel *Doc control II QA* *ztt 12 DEC 2024*
Name Title Signature/Date