

Product Number:	104570
Lot Number:	504307
Expiration:	6-Sep-2025
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104546	419104
Detector Reagent	104548	419105
SBG Reagent	104550	419106
Sample Diluent	104552	406619
Calibrator Diluent Reagent	104745	406620
Control Diluent Reagent	104554	406621
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

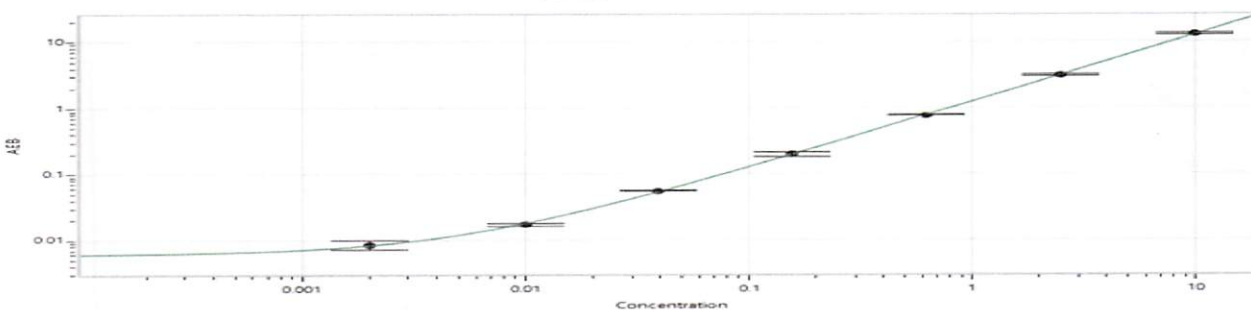
¹ Reagents are not Kit Lot Specific

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	
Control 1 Concentrate	406409	0.765	0.470 - 0.872
Control 2 Concentrate	406410	7.52	4.86 - 9.03

Calibrator Concentrate Reconstitution Volume (uL)	784
Control 1 Concentrate Reconstitution Volume (uL)	784
Control 2 Concentrate Reconstitution Volume (uL)	784

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

Calibration Curve



Review/Approval

Cody Carlson

QC Manager

Name

Title

[Signature] 01 Aug 2024
Signature/Date