

Product Number:	104570
Lot Number:	504180
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	406612
Detector Reagent	104548	406806
SBG Reagent	104550	406807
Sample Diluent	104552	406610
Calibrator Diluent	104745	406617
Control Diluent	104554	406618
Calibrator Concentrate	104566	406702
Control 1 Concentrate	104568	406703
Control 2 Concentrate	104569	406704
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

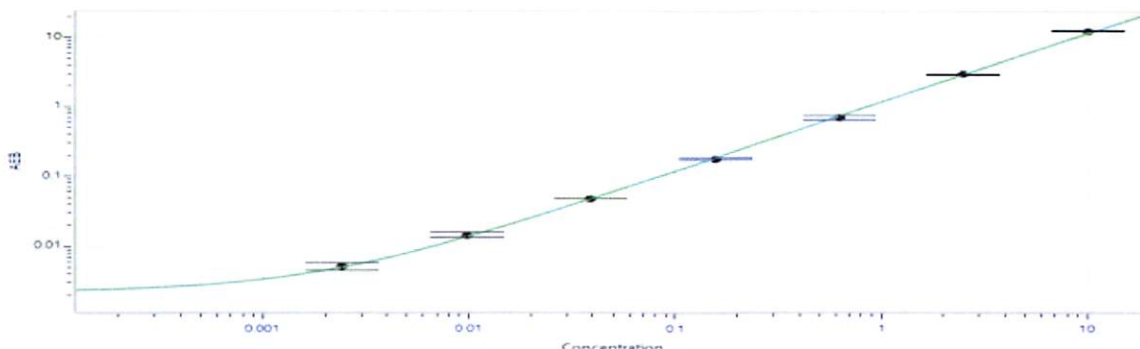
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	406703	0.663	0.466 - 0.866
Control 2 Concentrate	406704	7.18	5.06 - 9.39
Panel 1	405708	0.135	0.097 - 0.181
Panel 2	405708	9.37	7.06 - 13.1

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

Calibrator Concentrate Reconstitution Volume (uL)	649
Control 1 Concentrate Reconstitution Volume (uL)	649
Control 2 Concentrate Reconstitution Volume (uL)	649

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

Cody Carlson 25 APR 2024

Signature/Date