

Product Number:	104364
Lot Number:	504147
Expiration:	09-FEB-2026
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	104352	410803
Detector Reagent	104353	410702
SBG Reagent	104354	410703
Sample Diluent	104355	410704
Calibrator Diluent	104358	410614
Control Diluent	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

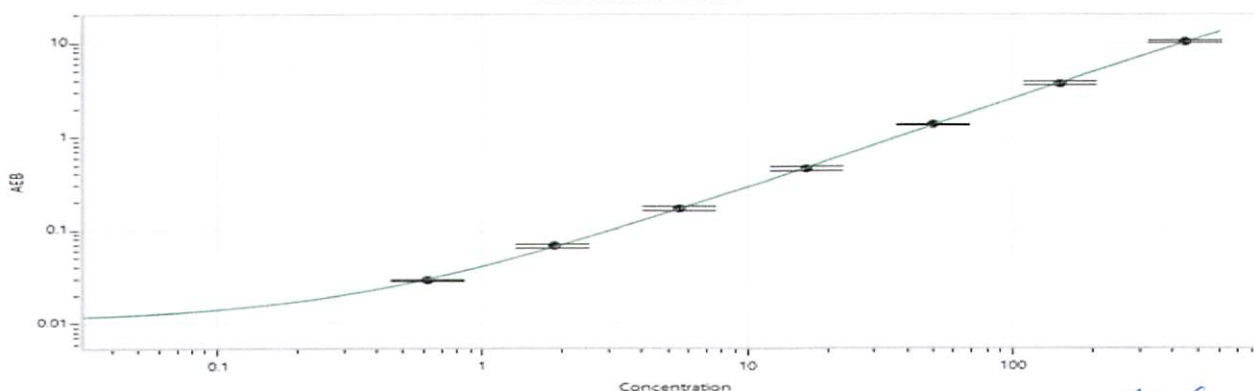
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	404520	4.27	3.48 - 6.19
Control 2	404521	352	265 - 524
Panel 1	409602	8.89	6.53 - 11.4
Panel 2	409602	4.83	4.61 - 6.92
Panel 3	409602	89.9	57.3 - 107
Panel 4	412201	386	188 - 442

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

Calibrator Concentrate Reconstitution Volume (uL)	1387
Control 1 Concentrate Reconstitution Volume (uL)	1387
Control 2 Concentrate Reconstitution Volume (uL)	1387

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

Calibration Curve



Review/Approval

Cody Carlson
Name Title

QC Manager

[Signature] 06 JUN 2024
Signature/Date