

Product Number:	104512
Lot Number:	504084
Expiration:	20-May-2025
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	104485	332407
Detector Reagent	104487	334513
SBG Reagent	104489	334107
Sample Diluent	104491	335602
Calibrator Diluent	104494	402505
Control Diluent	104495	402506
Calibrator Concentrate	104508	330613
Control 1 Concentrate	104510	330615
Control 2 Concentrate	104511	330616
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	330615	2.92	1.93 - 4.99
Control 2 Concentrate	330616	235	185 - 344
Panel 1	328515	43.5	27.6 - 51.3
Panel 2	328515	146	102 - 191
Panel 3	N/A	N/A	N/A
Panel 4	N/A	N/A	N/A

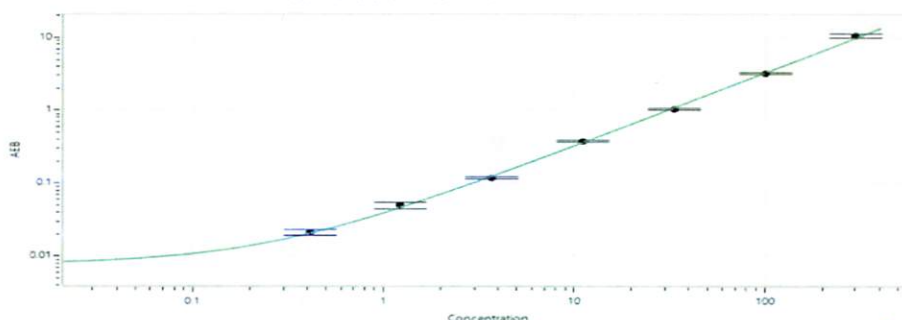
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	788
Control 1 Concentrate Reconstitution Volume (uL)	788
Control 2 Concentrate Reconstitution Volume (uL)	788

Calibrator Levels (pg/mL)

A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Calibration Curve



Review/Approval

Cody Carlson

QC Manager

Signature/Date

Name

Title

Signature/Date