

Product Number:	104618
Lot Number:	504098
Expiration:	18-Feb-2025
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Lot Number
Bead Reagent	104602	321411	N/A
Detector Reagent	104603	322208	N/A
SBG Reagent	104604	321316	N/A
Sample Diluent	104605	321317	N/A
Calibrator Diluent	104607	313811	313811A
Control Diluent	104608	313812	313812A
Calibrator Concentrate	104614	400901	N/A
Control 1 Concentrate	104616	400902	N/A
Control 2 Concentrate	104617	400903	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

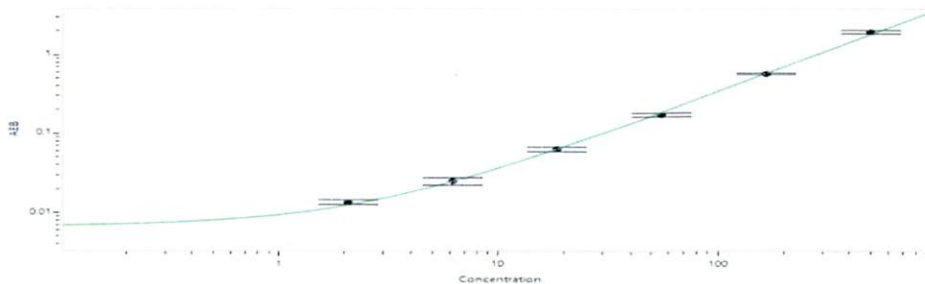
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	400902	62.1	49.8 - 89.8
Control 2 Concentrate	400903	1657	1468 - 2203
Panel 1	401906	14.6	11.7 - 19.7
Panel 2	401906	14.6	8.62 - 24.5
Panel 3	401906	729	604 - 905
Panel 4	N/A	N/A	N/A

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

Calibrator Concentrate Reconstitution Volume (uL)	1306
Control 1 Concentrate Reconstitution Volume (uL)	805
Control 2 Concentrate Reconstitution Volume (uL)	805

Calibrator Levels (pg/mL)	
A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500

Calibration Curve



Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 01 FEB 2024
Signature/Date