

Product Number:	Individual: 103769; Bulk: 103845	
Lot Number:	Individual: 504108; Bulk: 504109	
Expiration Date:	19-Jul-25	
Platform(s):	HD-X	
Component	Part Number	Lot Number
Bead Reagent	103758	404608
Detector Reagent	103760	404706
SBG Reagent	103762	404707
Sample Diluent	103763	404508
Calibrators	103766	405203
Control 1	103767	405108
Control 2	103768	405108
Control 3	103773	405108
RGP Reagent	103159	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 (ng/mL)	Mean Low Limit (ng/mL)	Mean High Limit (ng/mL)
Control 1	405108	324	245	369
Control 2	405108	2811	2120	3431
Control 3	405108	12225	10142	15212
	Lot	Result (ng/mL)	Mean Low Limit (ng/mL)	Mean High Limit (ng/mL)
Panel 1	319505	90.5	65.9	120
Panel 2	312101	64.3	49.2	125
Panel 3	312101	53.3	35.3	76.0
Panel 4	312101	34.2	28.4	46.7
Panel 5	312101	18.5	16.2	25.7
Panel 6	325503	8.20	7.09	10.6
Panel 7	312101	6.59	5.62	8.44
Panel 8	312101	4.84	4.03	6.04
Panel 9	312101	3.24	2.72	4.08
Panel 10	312101	2.28	1.91	2.86
Panel 11	312101	1.63	1.31	1.97
Panel 12	312101	1.11	0.847	1.27
Panel 13	312101	0.756	0.335	1.46
Panel 14	312101	0.618	0.499	0.823
Panel 15	312101	0.558	0.434	0.651
Panel 16	312101	0.381	0.318	0.494
Panel 17	312101	0.458	0.257	0.646
Panel 18	312101	0.020		
Panel 19	312101	0.032		
Panel 20	312101	0.072		

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

Review/Approval

Cody Carlson

QC Manager

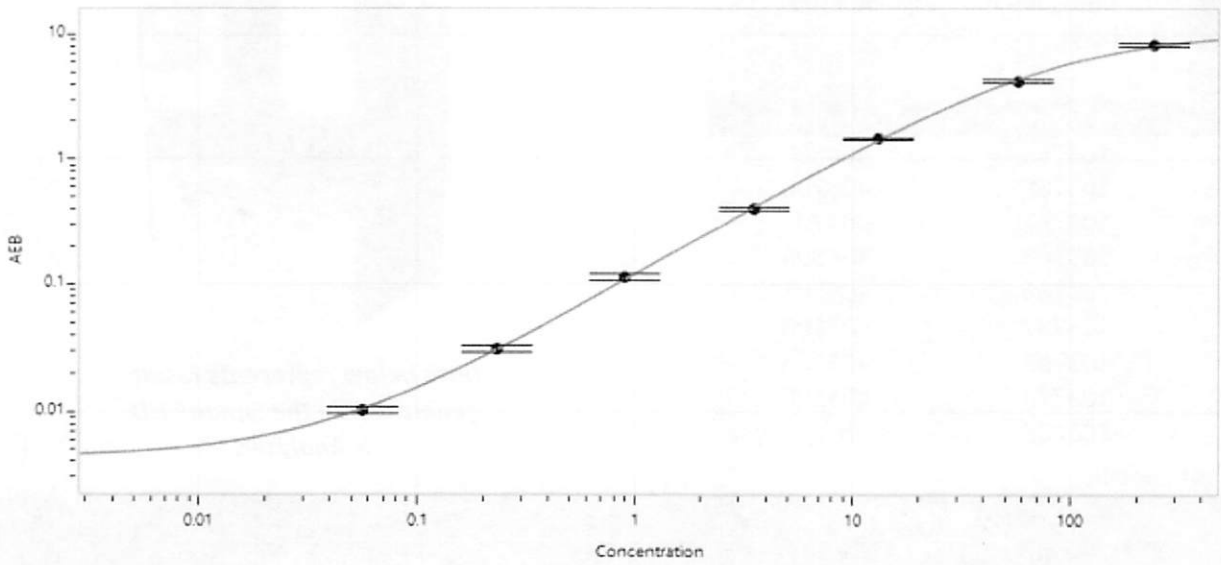
Name

Title

[Signature] 07MAR2024

Signature/Date

Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.056
C	0.235
D	0.892
E	3.50
F	13.1
G	56.9
H	242