

Product Number:	103769
Lot Number:	503657
Expiration:	22-Jun-2023
Platform(s):	HD-X
Product Number (Bulk):	103845
Lot Number (Bulk):	503661



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

Component	Part Number	Lot Number
Bead Reagent	103758	217313
Detector Reagent	103760	217605
SBG Reagent	103762	217606
Sample Diluent	103763	218005
Calibrators	103766	219205
Control 1	103767	232106
Control 2	103768	232106
Control 3	103773	232106
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				
Sample	Lot	Result (ng/mL)	Mean Low limit (ng/mL)	Mean High Limit (ng/mL)
Control 1	232106	350	229	561
Control 2	232106	3974	3067	4600
Control 3	232106	11606	9905	14858

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

Sample	Lot	Result (ng/mL)	Mean Low limit (ng/mL)	Mean High Limit (ng/mL)
Panel 1	121601	1.16	0.861	1.29
Panel 2	121601	2.22	1.60	2.39
Panel 3	121601	4.13	3.00	4.91
Panel 4	121601	7.54	5.74	9.12
Panel 5	121601	14.7	11.0	19.3
Panel 6	121601	28.3	20.2	35.7
Panel 7	121601	1.81	1.30	1.96
Panel 8	121601	6.72	3.74	8.16
Panel 9	121601	7.68	5.65	9.45
Panel 10	121601	18.5	12.6	21.6
Panel 11	121601	0.016	Reported Result	
Panel 12	121601	0.028	Reported Result	
Panel 13	121601	0.013	Reported Result	
Panel 14	121601	0.025	Reported Result	
Panel 15	121601	0.171	Reported Result	
Panel 16	121601	0.020	Reported Result	
Panel 17	121601	0.010	Reported Result	
Panel 18	121601	0.034	Reported Result	
Panel 19	121601	0.007	Reported Result	
Panel 20	121601	0.535	Reported Result	

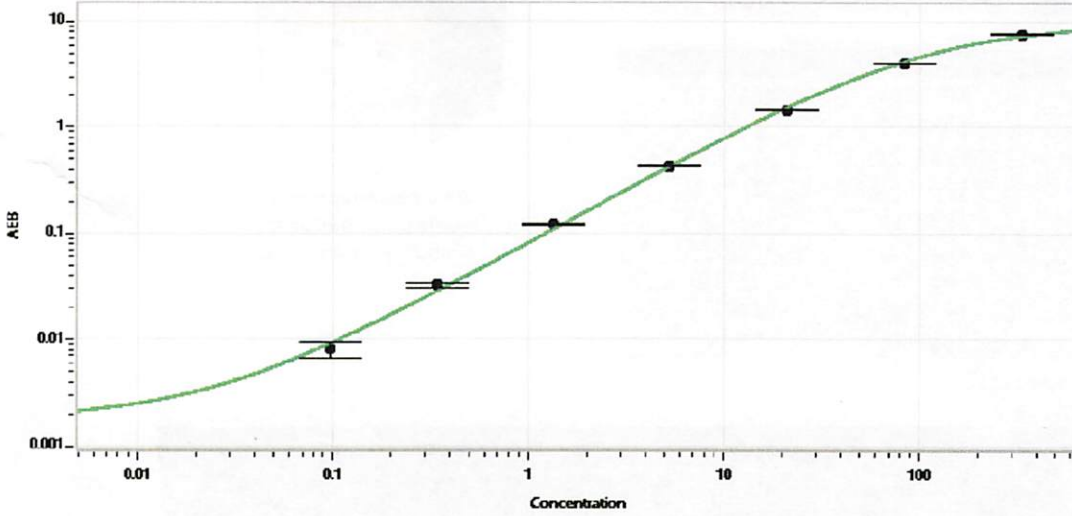
Review/Approval

Sonia Khiller
Name

ac Manager
Title

Sonia Khiller 15 DEC 2022
Signature/Date

Calibration Curve



Calibrator Levels
(ng/mL)

A	0.000
B	0.098
C	0.345
D	1.34
E	5.30
F	21.1
G	84.4
H	338