

Product Number:	103769
Lot Number:	503477
Expiration:	22-Jun-2023
Platform(s):	HD-X
Product Number (Bulk):	103845
Lot Number (Bulk):	503555



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	220201
Control 2	103768	220301
Control 3	103773	220301
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	220201	337	254	467
Control 2	220301	3685	2888	5065
Control 3	220301	16640	12270	23566

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

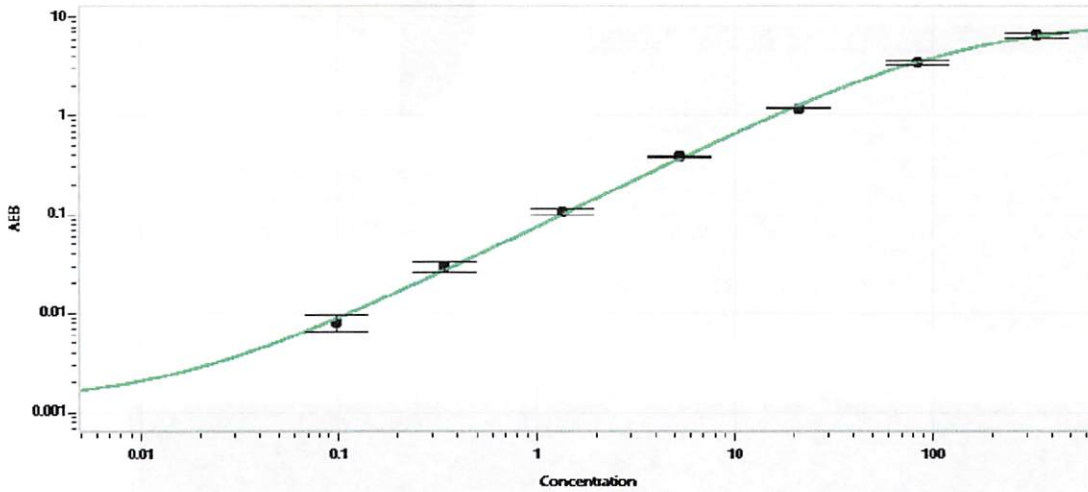
Sample	Lot	Result (ng/mL)	Mean Low limit (ng/mL)	Mean High Limit (ng/mL)
Panel 1	121601	1.14	0.861	1.29
Panel 2	121601	1.84	1.60	2.39
Panel 3	121601	4.01	3.00	4.91
Panel 4	121601	7.95	5.74	9.12
Panel 5	121601	15.4	11.0	19.3
Panel 6	121601	31.7	20.2	35.7
Panel 7	121601	1.80	1.30	1.96
Panel 8	121601	6.73	3.74	8.16
Panel 9	121601	8.70	5.65	9.45
Panel 10	121601	19.6	12.6	21.6
Panel 11	121601	0.015		Reported Result
Panel 12	121601	0.022		Reported Result
Panel 13	121601	0.016		Reported Result
Panel 14	121601	0.015		Reported Result
Panel 15	121601	0.163		Reported Result
Panel 16	121601	0.020		Reported Result
Panel 17	121601	0.017		Reported Result
Panel 18	121601	0.009		Reported Result
Panel 19	121601	0.019		Reported Result
Panel 20	121601	0.575		Reported Result

Review/Approval

Sonia Khillan QC Manager
Name Title

Sonia Khillan / 01/NOV2022
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