

Product Number:	103769
Lot Number:	503444
Expiration:	27-Feb-2023
Platform(s):	HD-X
Product Number (Bulk):	103845
Lot Number (Bulk):	N/A



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

Component	Part Number	Lot Number
Bead Reagent	103758	211201
Detector Reagent	103760	214706
SBG Reagent	103762	215401
Sample Diluent	103763	212805
Calibrators	103766	213009
Control 1	103767	214104
Control 2	103768	214104
Control 3	103773	214104
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	214104	462	241	524
Control 2	214104	4336	2902	4353
Control 3	214104	21351	14506	21759

* 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	1.18	0.846	1.28
Panel 2	121601	2.25	1.60	2.40
Panel 3	121601	4.15	2.91	4.85
Panel 4	121601	7.58	5.62	8.80
Panel 5	121601	14.9	10.3	19.9
Panel 6	121601	27.0	20.6	32.8
Panel 7	121601	1.75	1.28	1.92
Panel 8	121601	6.35	3.84	7.36
Panel 9	121601	8.06	5.77	8.80
Panel 10	121601	19.2	12.9	19.8
Panel 11	121601	0.018	Reported Result	
Panel 12	121601	0.022	Reported Result	
Panel 13	121601	0.024	Reported Result	
Panel 14	121601	0.014	Reported Result	
Panel 15	121601	0.197	Reported Result	
Panel 16	121601	0.033	Reported Result	
Panel 17	121601	0.028	Reported Result	
Panel 18	121601	0.044	Reported Result	
Panel 19	121601	0.038	Reported Result	
Panel 20	121601	0.575	Reported Result	

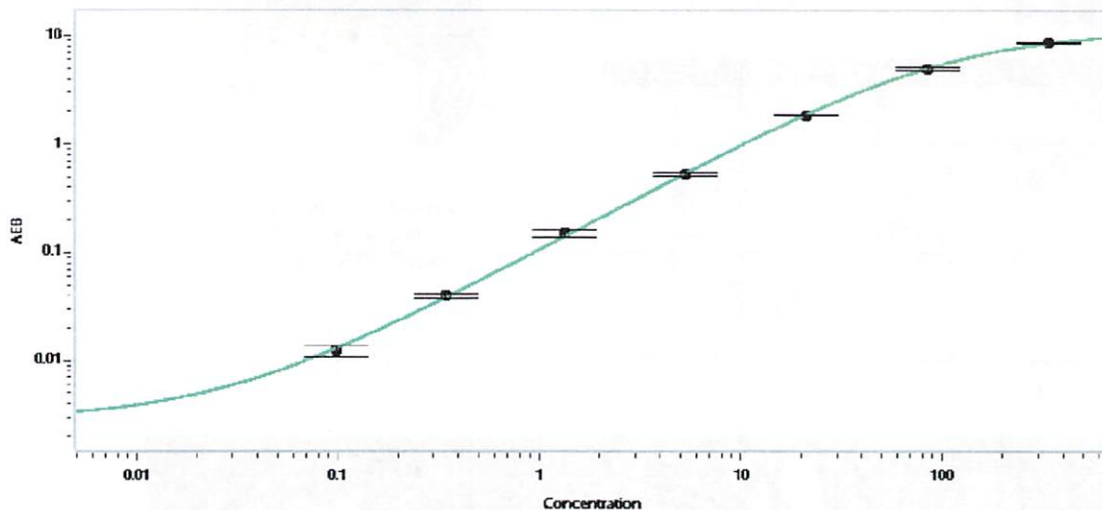
Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 14 JUL 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