

Product Number:	103769
Lot Number:	502764
Expiration:	6-Oct-2021
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	028007
Detector Reagent	103760	102901
SBG Reagent	103762	106819
Sample Diluent	103763	103414
Calibrators	103766	103503
Control 1	103767	103504
Control 2	103768	103504
Control 3	103773	103504
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	103504	419	194	456
Control 2	103504	2346	2220	3567
Control 3	103504	14193	11921	17960

* 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	020901	17.4	12.7	23.1
Panel 2	020901	4.55	3.88	5.82
Panel 3	020901	20.5	15.0	22.5
Panel 4	020901	15.1	13.7	20.9
Panel 5	020901	21.5	16.7	30.7
Panel 6	020901	19.3	13.9	20.8
Panel 7	020901	3.89	2.83	4.55
Panel 8	020901	1.28	1.03	1.61
Panel 9	020901	3.96	2.90	5.02
Panel 10	020901	78.6	40.3	115
Panel 11	020901	0.309	Reported Result	
Panel 12	020901	0.065	Reported Result	
Panel 13	020901	0.090	Reported Result	
Panel 14	020901	0.018	Reported Result	
Panel 15	020901	0.029	Reported Result	
Panel 16	020901	0.054	Reported Result	
Panel 17	031107	0.051	Reported Result	
Panel 18	031109	0.017	Reported Result	
Panel 19	106401	0.040	Reported Result	
Panel 20	106402	0.006	Reported Result	

Review/Approval

Junpil Park QC Supervisor

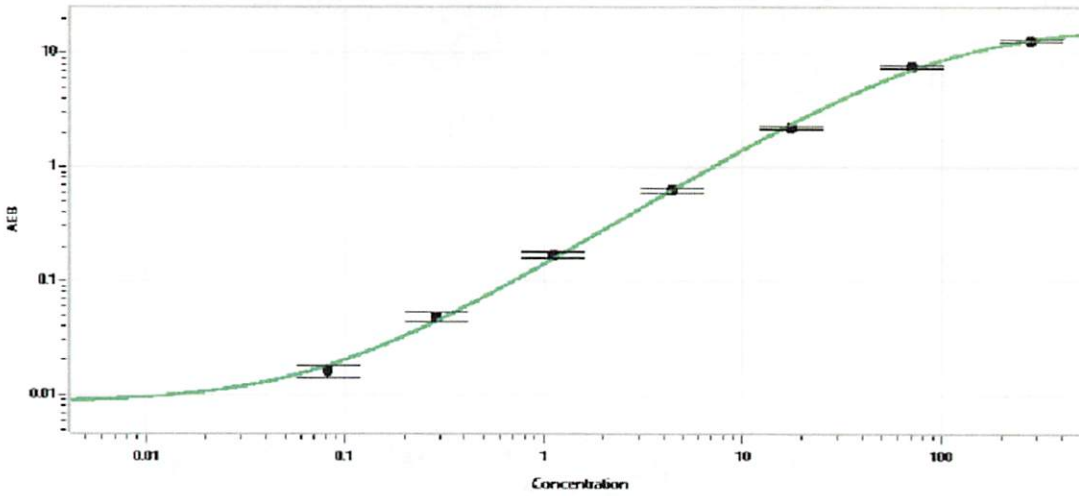
Name

Title

31 Mar 2021

Signature/Date

Calibration Curve



**Calibrator Levels
(ng/mL)**

A	0.000
B	0.049
C	0.196
D	0.783
E	3.13
F	12.5
G	50.1
H	224