

Product Number:	103769
Lot Number:	503065
Expiration:	11-Jun-2022
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	125701
Detector Reagent	103760	126612
SBG Reagent	103762	125404
Sample Diluent	103763	125502
Calibrators	103766	125803
Control 1	103767	125804
Control 2	103768	125804
Control 3	103773	125804
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	125804	397	217	540
Control 2	125804	3653	2879	4318
Control 3	125804	17392	10800	23283

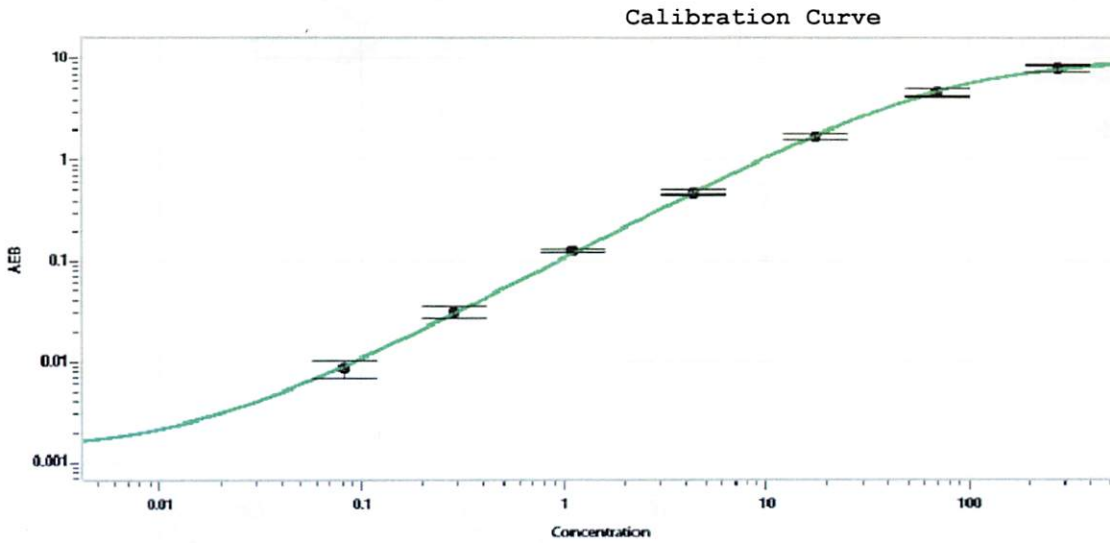
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Sample	Lot	Result (ng/mL)	Mean Low limit (ng/mL)	Mean High Limit (ng/mL)
Panel 1	111710	1.10	0.786	1.30
Panel 2	111710	1.94	1.46	2.31
Panel 3	111710	3.73	2.92	4.47
Panel 4	111710	7.21	5.57	8.35
Panel 5	111710	15.8	9.63	18.1
Panel 6	111710	28.9	18.2	35.2
Panel 8	111710	7.11	4.81	8.06
Panel 9	111710	15.0	6.84	16.2
Panel 10	111710	56.8	23.0	67.1
Panel 11	111710	0.249	Reported Result	
Panel 12	111710	0.047	Reported Result	
Panel 13	111710	0.031	Reported Result	
Panel 14	111710	0.033	Reported Result	
Panel 15	111710	0.045	Reported Result	
Panel 16	111710	0.035	Reported Result	
Panel 17	111710	0.054	Reported Result	
Panel 18	111710	0.033	Reported Result	
Panel 19	111710	0.018	Reported Result	
Panel 20	111710	0.539	Reported Result	

Review/Approval

Junpil Park    QC Supervisor  
Name                      Title

 29 Sep 2021  
Signature/Date



Calibrator Levels (ng/mL)	
A	0.000
B	0.082
C	0.287
D	1.11
E	4.39
F	17.5
G	69.9
H	280