

Product Number:	103769
Lot Number:	502935
Expiration:	8-May-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	112805
Detector Reagent	103760	113414
SBG Reagent	103762	112806
Sample Diluent	103763	112807
Calibrators	103766	107607
Control 1	103767	107804
Control 2	103768	107804
Control 3	103773	107804
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	107804	525	234	609
Control 2	107804	3844	2937	4406
Control 3	107804	16442	11941	20550

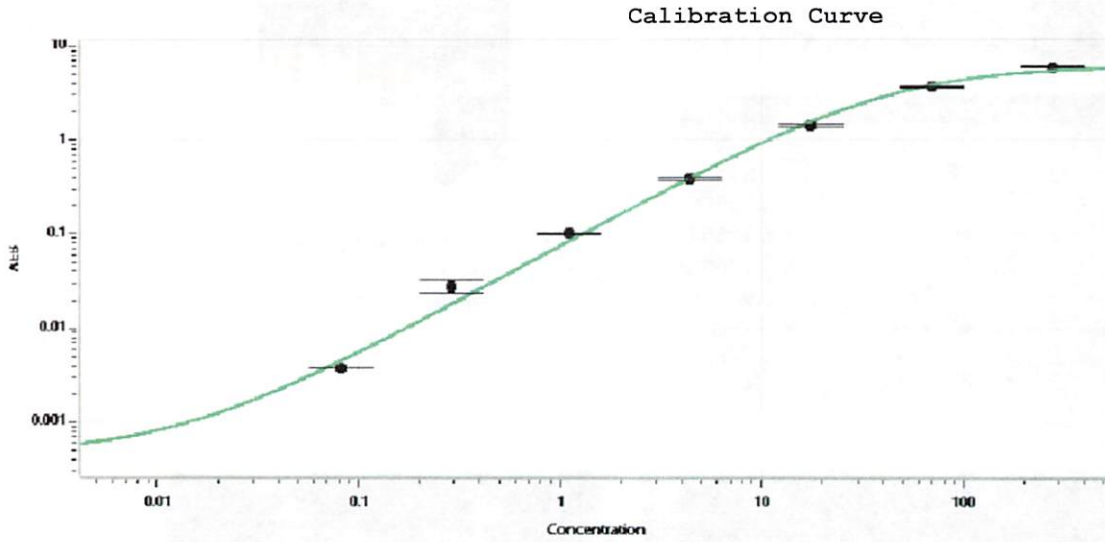
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Sample	Lot	Result (ng/mL)	Mean Low limit (ng/mL)	Mean High Limit (ng/mL)
Panel 1	111710	1.17	0.707	1.26
Panel 2	111710	2.00	1.39	2.08
Panel 3	111710	3.79	2.75	4.13
Panel 4	111710	6.55	5.27	7.90
Panel 5	111710	12.4	9.65	17.0
Panel 6	111710	23.3	17.4	35.3
Panel 7	111710	12.8	7.68	23.1
Panel 8	111710	6.44	4.87	7.30
Panel 9	111710	10.1	8.35	12.5
Panel 10	111710	40.2	20.4	66.3
Panel 11	111710	0.373	Reported Result	
Panel 12	111710	0.118	Reported Result	
Panel 13	111710	0.106	Reported Result	
Panel 14	111710	0.112	Reported Result	
Panel 15	111710	0.145	Reported Result	
Panel 16	111710	0.140	Reported Result	
Panel 17	111710	0.202	Reported Result	
Panel 18	111710	0.091	Reported Result	
Panel 19	111710	0.056	Reported Result	
Panel 20	111710	0.668	Reported Result	

Review/Approval

Junpil Park QC Supervisor
Name Title

 16 Jun 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