

Product Number:	103806
Lot Number:	502962
Expiration:	23-May-2022
Bulk Kit	
Product Number:	103872
Lot Number:	502963
Platform(s):	HD-X



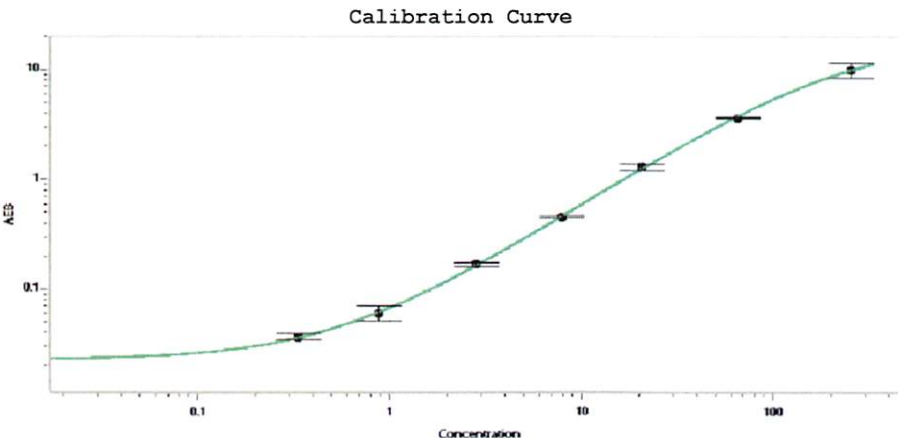
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103794	114406
Detector Reagent	103796	114309
SBG Reagent	103798	114408
Sample Diluent	103800	114410
Calibrators	103803	114519
Control 1	103804	114520
Control 2	103805	114520
RGP Reagent	103159	N/A ¹


¹ RGP is not Kit Lot Specific

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	114520	7.96	4.04-11.5
Control 2	114520	158	131-207
Panel 1	021305	2.74	2.55-4.27
Panel 2	021111	62.0	42.2-90.4
Panel 3	021111	123	89.1-202

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



Calibrator Levels (pg/mL)	
A	0.000
B	0.338
C	0.885
D	2.85
E	7.88
F	20.5
G	65.1
H	252

Review/Approval Junpil Park QC Supervisor  01 Jul 2021
Name Title Signature/Date