

Product Number:	103806
Lot Number:	502744
Expiration:	11-Jan-2022
	Bulk Kit
Product Number:	103872
Lot Number:	502745
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103794	104203
Detector Reagent	103796	104011
SBG Reagent	103798	105517
Sample Diluent	103800	104018
Calibrators	103803	104703
Control 1	103804	104705
Control 2	103805	104705
RGP Reagent	103159	N/A ¹

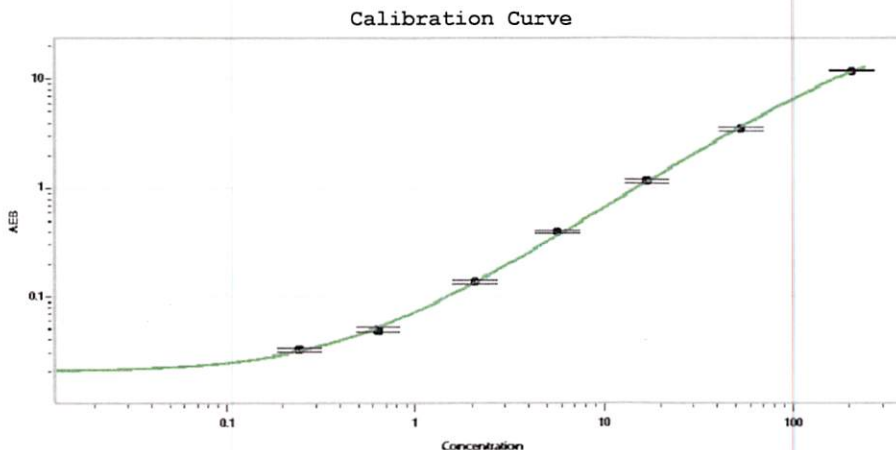
¹ RGP is not Kit Lot Specific



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

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	104705	6.67	5.04-7.56
Control 2	104705	136	110-172
Panel 1	021305	3.58	2.55-4.27
Panel 2	021111	51.9	42.2-90.4
Panel 3	021111	132	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	0.245
C	0.641
D	2.07
E	5.71
F	16.7
G	53.2
H	206

Review/Approval

Junpil Park QC Supervisor
Name Title

Signature/Date

[Signature] 16 Mar 2021