

Product Number:	103806
Lot Number:	502864
Expiration:	7-Apr-2022
	Bulk Kit
Product Number:	103872
Lot Number:	502865
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103794	110107
Detector Reagent	103796	110110
SBG Reagent	103798	110112
Sample Diluent	103800	110114
Calibrators	103803	109607
Control 1	103804	109608
Control 2	103805	109608
RGP Reagent	103159	N/A ¹

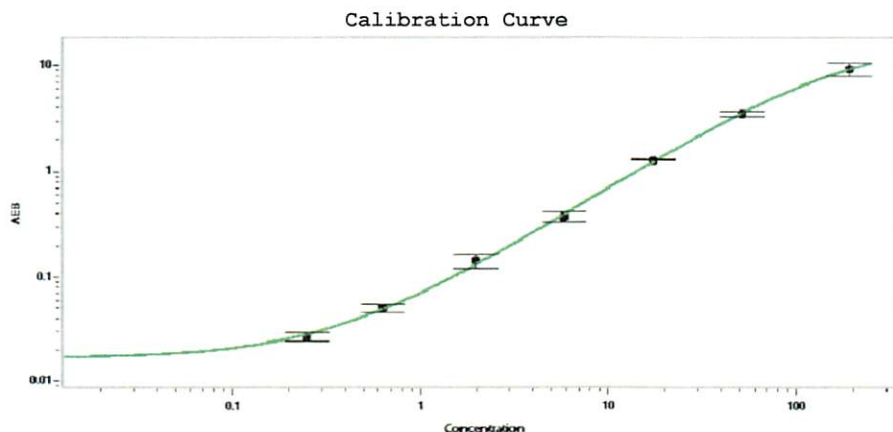
¹ RGP is not Kit Lot Specific



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

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	109608	6.48	4.10-6.81
Control 2	109608	118	96.0-166
Panel 1	021305	2.78	2.55-4.27
Panel 2	021111	47.2	42.2-90.4
Panel 3	021111	105	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	0.250
C	0.632
D	1.97
E	5.79
F	17.3
G	51.8
H	190

Review/Approval

Junpil Park QC Supervisor
Name Title

Signature/Date

[Signature] 29 Apr 2021