

Product Number:	103806
Lot Number:	502708
Expiration:	1-Dec-2021
Bulk Kit	
Product Number:	103872
Lot Number:	502710
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103794	102514
Detector Reagent	103796	102518
SBG Reagent	103798	102607
Sample Diluent	103800	102524
Calibrators	103803	102601
Control 1	103804	102602
Control 2	103805	102602
RGP Reagent	103159	N/A ¹

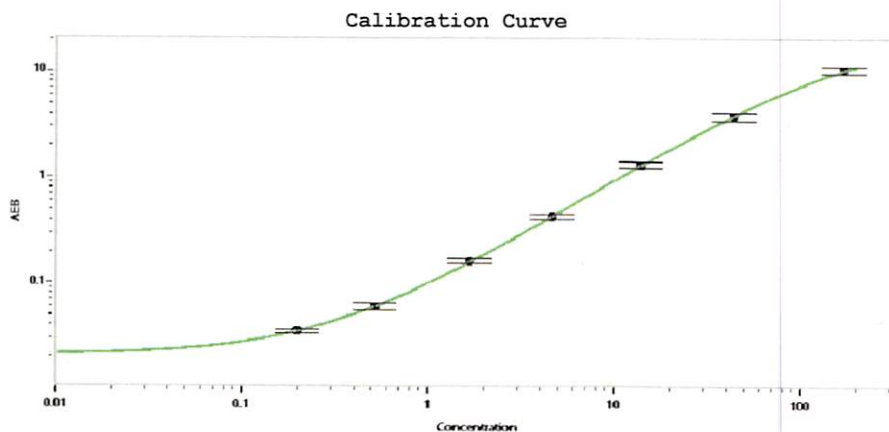
¹ RGP is not Kit Lot Specific



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

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	102602	5.11	4.25-6.38
Control 2	102602	103	89.0-146
Panel 1	021305	3.24	2.55-4.27
Panel 2	021111	59.0	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	0.199
C	0.521
D	1.68
E	4.64
F	13.8
G	43.9
H	170

Review/Approval Junpil Park QC Supervisor
Name Title

 12 Feb 2021
Signature/Date