

Product Number:	103806
Lot Number:	502679
Expiration:	15-Dec-2021
	Bulk Kit
Product Number:	103872
Lot Number:	502680
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103794	036503
Detector Reagent	103796	036502
SBG Reagent	103798	101212
Sample Diluent	103800	036501
Calibrators	103803	036401
Control 1	103804	036402
Control 2	103805	036402
RGP Reagent	103159	N/A ¹

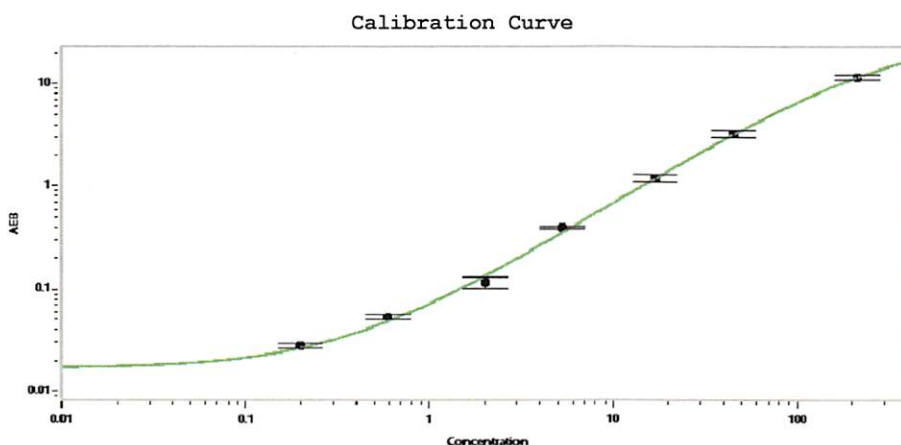
¹ RGP is not Kit Lot Specific



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

Release Materials*			
Sample	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	036402	6.39	4.32-7.50
Control 2	036402	141	102-189
Panel 1	021305	3.61	2.56-4.35
Panel 2	021111	69.2	43.1-93.0
Panel 3	021111	152	97.0-205

* 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	0.197
C	0.595
D	2.01
E	5.29
F	17.0
G	45.1
H	213

Review/Approval

Junpil Park QC Supervisor
Name Title

Signature/Date

25 Jan 2021