

Product Number:	101552
Lot Number:	503794
Expiration:	13-Jan-2024
Platform(s):	HD-1/HD-x & SR-X

Component	Part Number	Lot Number
Bead Reagent	101554	304007
Detector Reagent	101553	304106
SBG Reagent	101555	304408
Sample Diluent	101556	304409
Calibrators	101559	304601
Control 1	101560	304602
Control 2	101561	304602
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

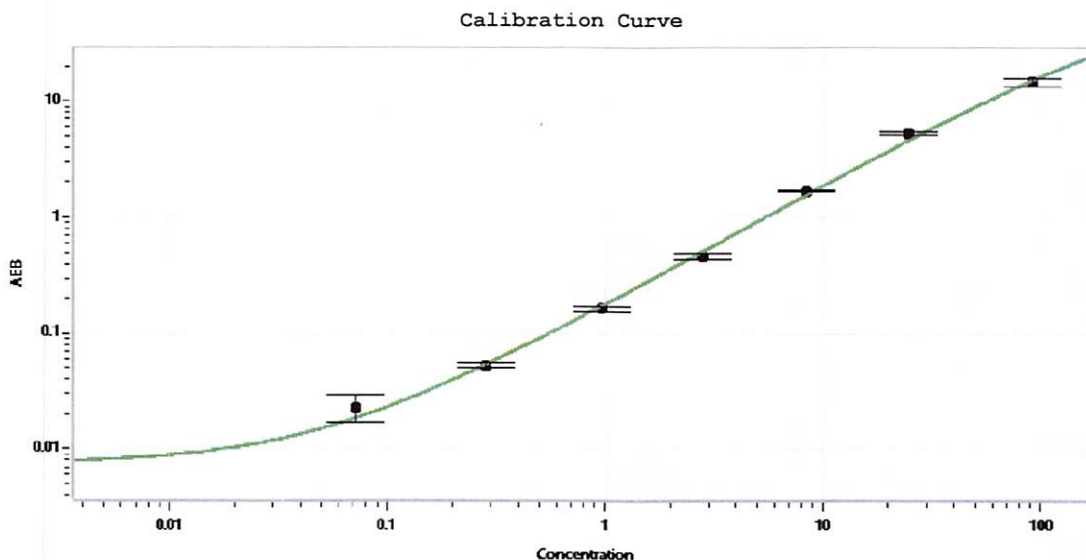
¹ RGP is not Kit Lot Specific



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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	304602	1.24	0.882-1.52
Control 2	304602	62.1	49.3-74.0
Panel 1	232607	1.14	0.576-1.41
Panel 2	232607	1.004	0.544-1.42
Panel 3	234908	22.5	16.7-25.0
Panel 4	234908	31.4	25.6-38.4

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



Calibrator Levels (pg/mL)	Value
A	0.000
B	0.119
C	0.335
D	1.10
E	3.13
F	10.1
G	30.6
H	88.0

Review/Approval: ROBERT SOMOGYI QC, MANAGER [Signature] 21 APR 2013
 Name Title Signature/Date