

Product Number:	101552
Lot Number:	502310
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-x & SR-X



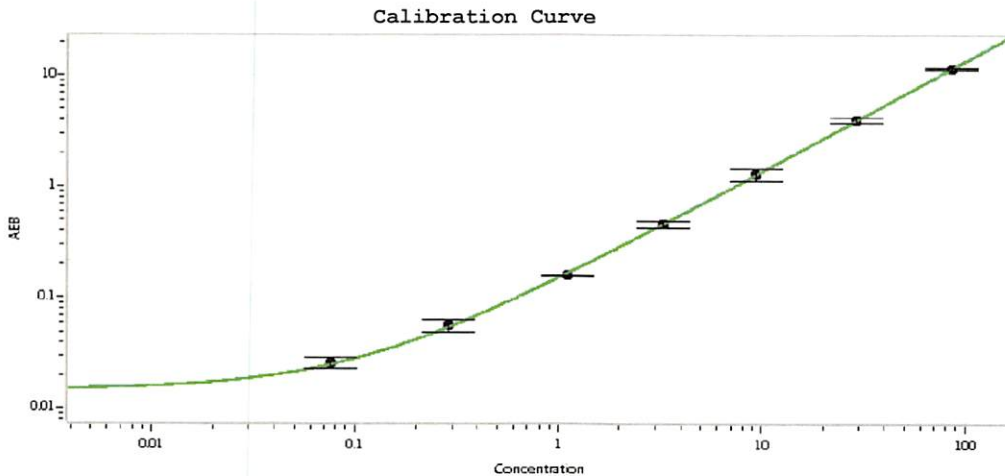
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101554	008003
Detector Reagent	101553	008004
SBG Reagent	101555	008304
Sample Diluent	101556	007809
Calibrators	101559	008601
Controls	101560-101561	008602
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008602	1.10	0.464-1.57
Control 2	008602	61.8	46.4-69.6
Panel 1	916901	0.777	0.089-1.59
Panel 2	916902	1.32	0.538-2.13
Panel 3	916903	29.6	15.5-33.4
Panel 4	916904	97.0	63.0-102

* 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	0.076
C	0.289
D	1.10
E	3.24
F	9.33
G	28.7
H	85.5

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 06 Apr 2020
Signature/Date