

Product Number:	101552
Lot Number:	501903
Expiration:	13-Jun-2020
Platform(s):	HD-1/HD-x & SR-X



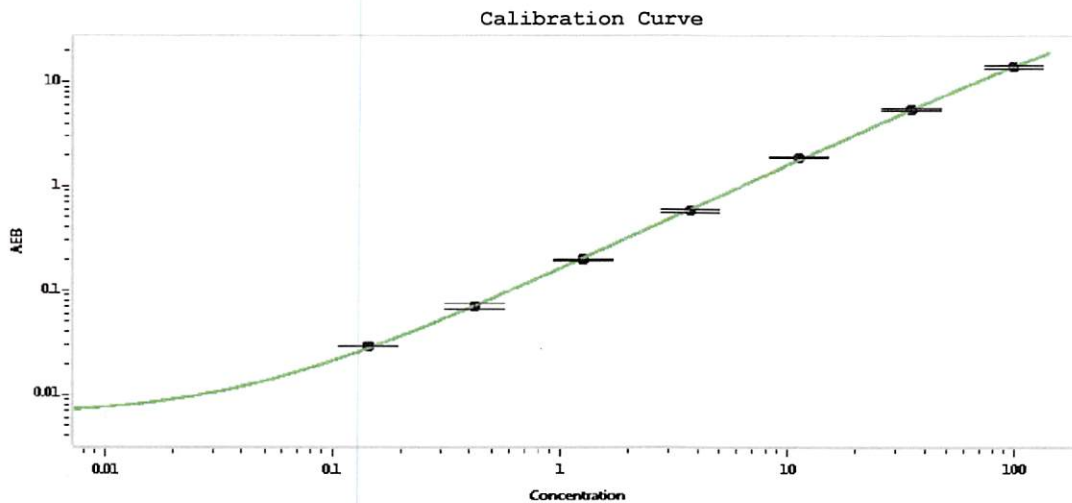
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101554	918210
Detector Reagent	101553	918206
SBG Reagent	101555	918205
Sample Diluent	101556	918207
Calibrators	101559	918404
Controls	101560-101561	918405
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	918405	0.914	0.718-1.82
Control 2	918405	58.3	47.6-71.5
Panel 1	916901	0.992	0.691-1.40
Panel 2	916902	1.55	1.22-2.00
Panel 3	916903	24.1	17.1-28.8
Panel 4	916904	83.3	62.7-94.0

* 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	0.088
C	0.315
D	1.01
E	3.00
F	9.55
G	27.9
H	91.1

Review/Approval

Patrizia Stadler Assoc. Director, QC *P. Stadler* 09 OCT 2019
 Name Title Signature/Date