

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



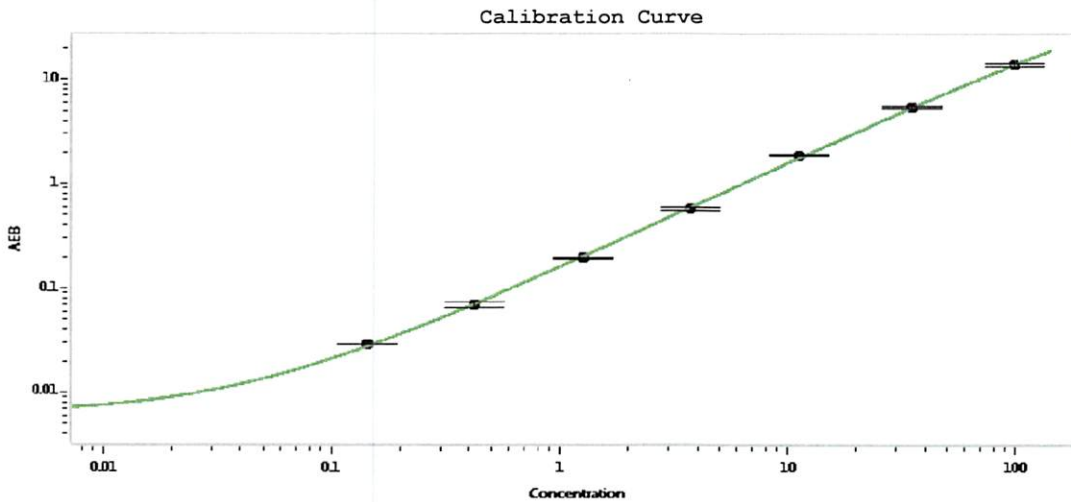
Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	1.26	0.897-1.74
Control 2	918405	68.7	48.8-73.2
Panel 1	916901	1.37	0.619-1.52
Panel 2	916902	2.00	1.17-2.09
Panel 3	916903	28.3	17.0-29.8
Panel 4	916904	93.8	64.0-96.0

* 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	0.143
C	0.422
D	1.27
E	3.71
F	11.2
G	35.1
H	98.6

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P Stadler 25 Jul 2019
Signature/Date