

Product Number:	100860
Lot Number:	501929
Expiration:	6-Mar-2020
Platform(s):	HD-1/ HD-X



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

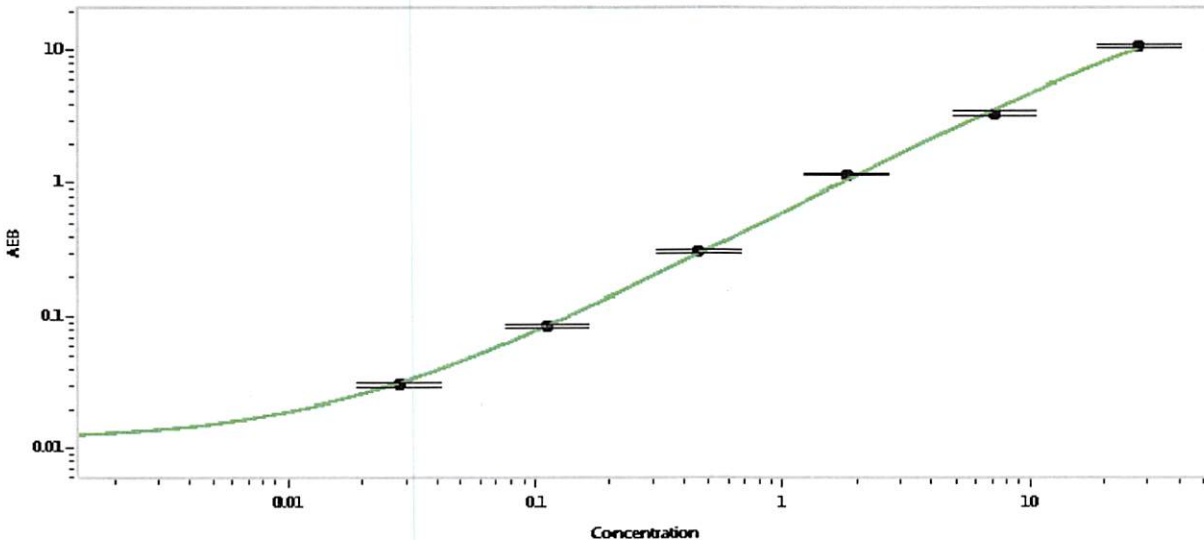
Component	Part Number	Lot Number
Bead Reagent	100865	908706
Detector Reagent	100862	906508
SBG Reagent	100868	906507
Sample Diluent	100870	906506
Calibrators	101111	907103
Controls	100875 & 100876	910602
RGP Reagent	103159	N/A1

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910602	6.29	3.53-10.6
Control 2	910602	0.117	0.061-0.182
Panel 1	908501	1.68	1.30-1.94
Panel 2	908502	9.37	7.54-12.2
Panel 3	908503	39.0	23.6-46.4

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.028
C	0.112
D	0.459
E	1.82
F	7.20
G	27.3

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	18 Sep 2019
	Name	Title	Signature/Date	