

Product Number:	100860
Lot Number:	501799
Expiration:	16-Apr-2020
Platform(s):	HD-1



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

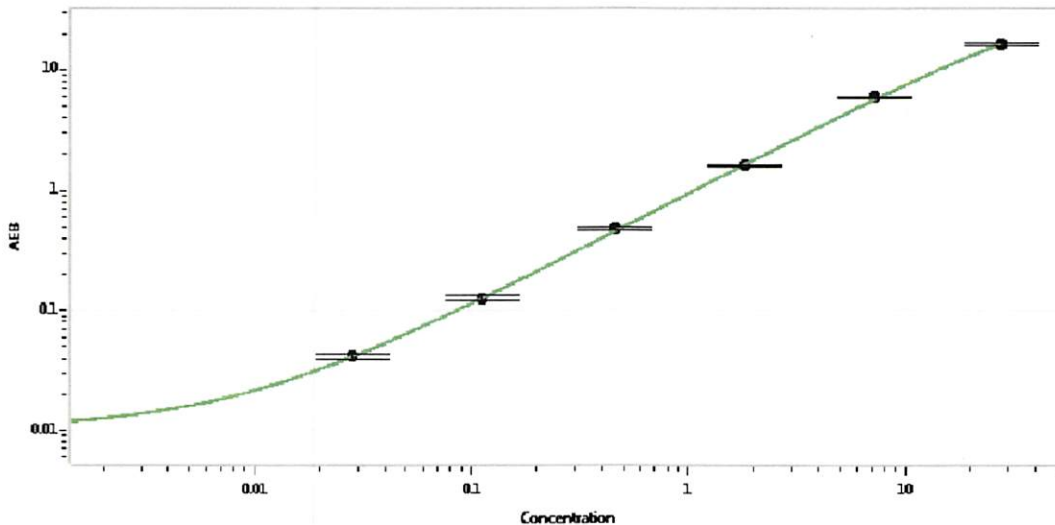
Component	Part Number	Lot Number
Bead Reagent	100865	911410
Detector Reagent	100862	911506
SBG Reagent	100868	911311
Sample Diluent	100870	911310
Calibrators	101111	907103
Controls	100875-100876	910602
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910602	6.11	3.46-10.4
Control 2	910602	0.110	0.060-0.180
Panel 1	908501	1.40	1.27-1.91
Panel 2	908502	9.65	7.66-12.0
Panel 3	908503	34.1	24.2-45.5

* 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	27.30

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	01MAY 2019
	Name	Title	Signature/Date	