

Product Number:	100860
Lot Number:	502016
Expiration:	6-Jun-2020
Platform(s):	HD-1/ HD-X



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

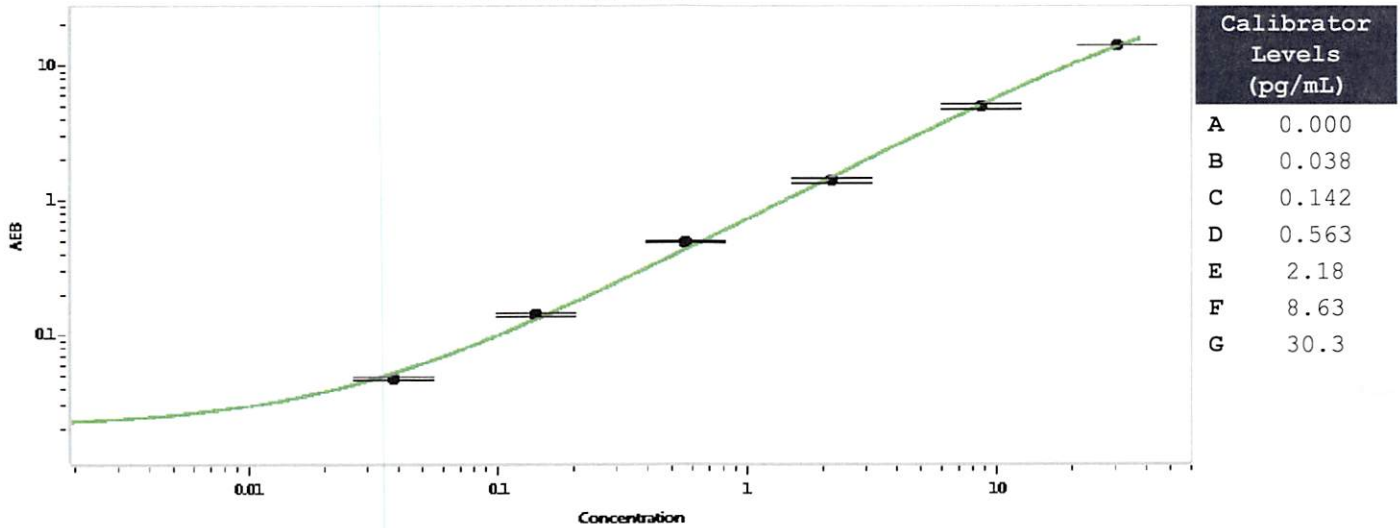
Component	Part Number	Lot Number
Bead Reagent	100865	924608
Detector Reagent	100862	919602
SBG Reagent	100868	920706
Sample Diluent	100870	915705
Calibrators	101111	923904
Controls	100875-100876	923905
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	923905	6.27	2.97-8.90
Control 2	923905	0.113	0.052-0.157
Panel 1	908501	1.45	1.17-1.91
Panel 2	908502	8.50	7.29-12.0
Panel 3	908503	30.9	23.8-44.9

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

Calibration Curve



Review/Approval: Patrizia Stadler Assoc. Director, QC *P. Stodola* 17 Sep 2019
Name Title Signature/Date