

Product Number:	100860
Lot Number:	502084
Expiration:	3-Sep-2020
Platform(s):	HD-1/ HD-X



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

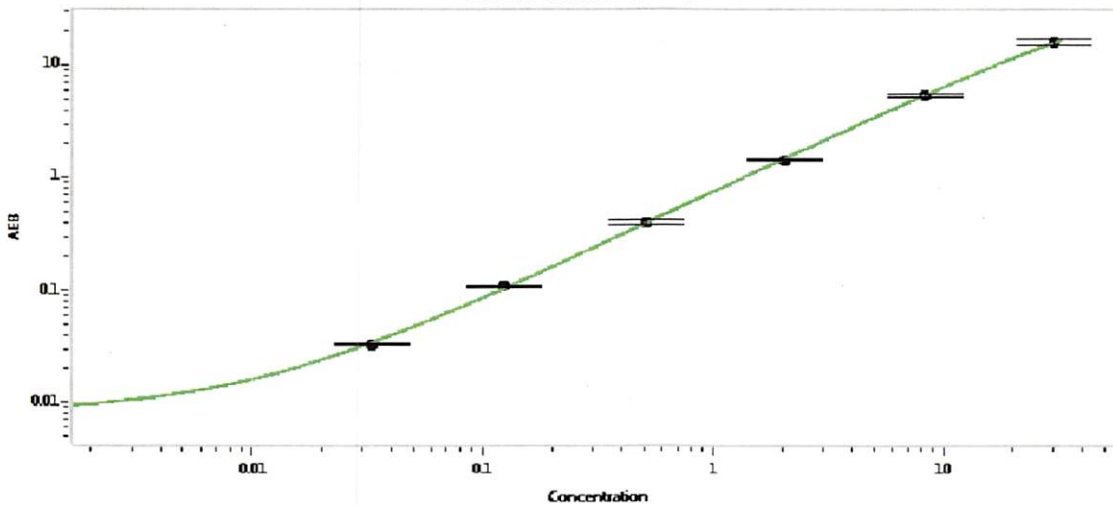
Component	Part Number	Lot Number
Bead Reagent	100865	926805
Detector Reagent	100862	929001
SBG Reagent	100868	924607
Sample Diluent	100870	924610
Calibrators	101111	926101
Controls	100875 & 100876	926102
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	926102	5.36	4.42-6.63
Control 2	926102	0.080	0.069-0.103
Panel 1	908501	1.09	0.978-1.95
Panel 2	908502	7.61	6.73-11.8
Panel 3	908503	24.9	20.5-44.6

\* 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	0.033
C	0.124
D	0.512
E	2.03
F	8.20
G	30.1

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p. stadler</i> 14 Nov 2019
	Name	Title	Signature/Date