

Product Number:	100860
Lot Number:	502346
Expiration:	30-Jan-2021
Platform(s):	HD-1/ HD-X



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

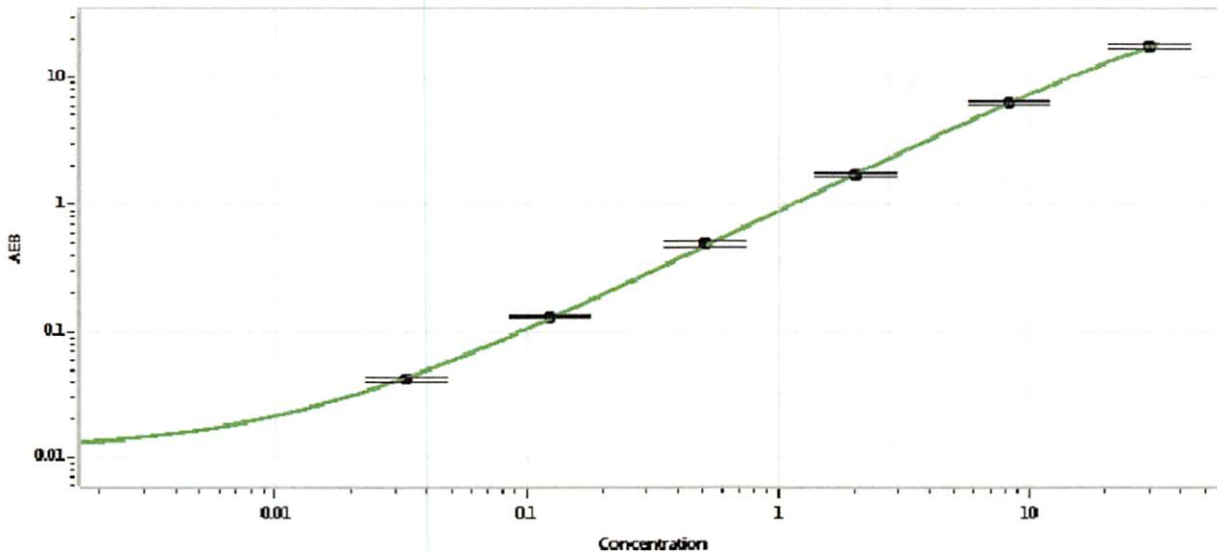
Component	Part Number	Lot Number
Bead Reagent	100865	010712
Detector Reagent	100862	007206
SBG Reagent	100868	011110
Sample Diluent	100870	003004
Calibrators	101111	926101
Controls	100875 & 100876	926102
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	926102	6.00	4.48-6.71
Control 2	926102	0.091	0.068-0.109
Panel 1	908501	0.926	0.631-2.10
Panel 2	908502	6.80	4.38-12.5
Panel 3	908503	22.9	12.1-47.2

* 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	Shivani Goel Name	QC Manager Title	Shivani Goel 29 Apr 2020 Signature/Date
-----------------	----------------------	---------------------	--------------------------------------------