

Product Number:	101235
Lot Number:	502400
Expiration:	7-Oct-2020
Platform(s):	HD-1/HD-X



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

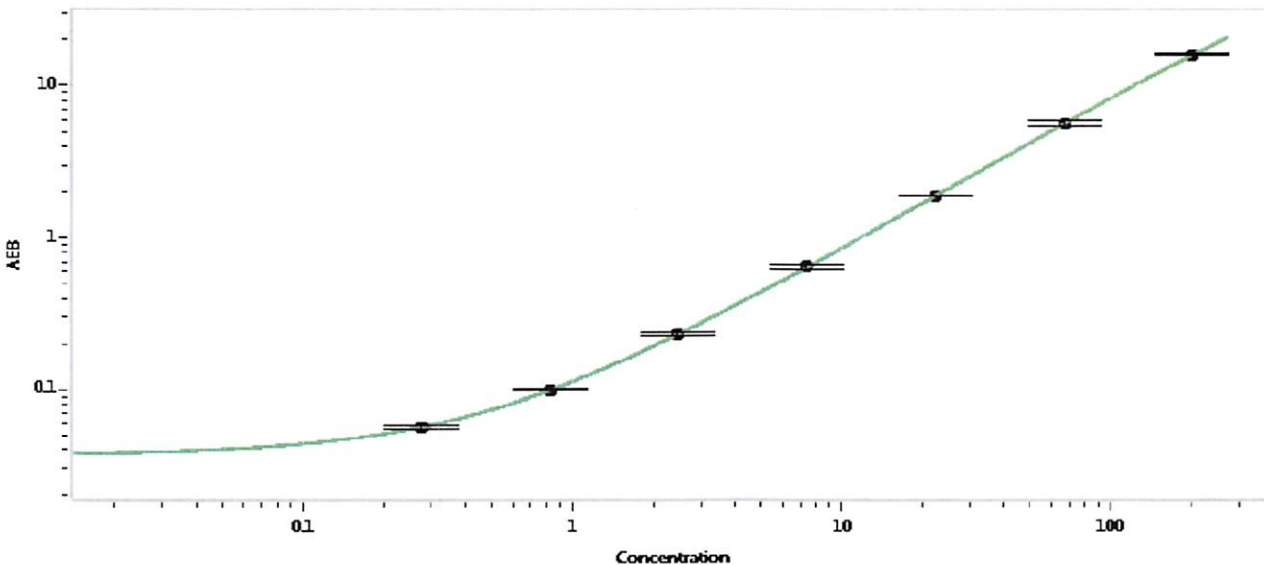
Component	Part Number	Lot Number
Bead Reagent	101222	930905
Detector Reagent	101224	930315
SBG Reagent	101234	931005
Sample Diluent	101232	930209
Calibrator Diluent	101228	928022
Calibrator Concentrate	101226	931903
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	62432	16.6	13.2-19.8
Control 2	62432	460	396-593
Panel 1	MF-081516-01	8.60	5.69-10.1
Panel 2	MF-081516-02	87.8	52.3-90.4
Panel 3	MF-081516-03	3.03	1.54-4.72

* 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	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shwam Goel 02 Jun 2020
Signature/Date