

Product Number:	101664
Lot Number:	502468
Expiration:	22-Jan-2021
Platform(s):	HD-1/HD-X



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

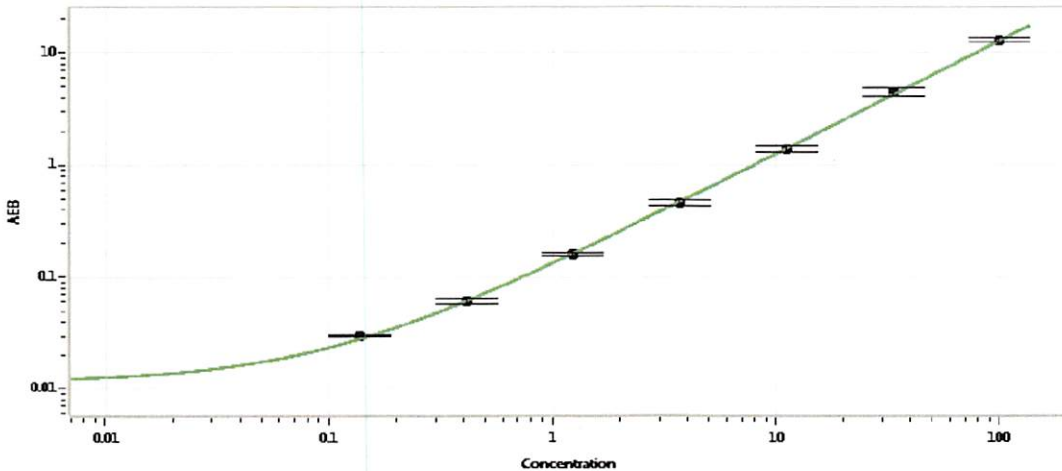
Component	Part Number	Lot Number
Bead Reagent	101666	002215
Detector Reagent	101665	002203
SBG Reagent	101667	019009
Sample Diluent	101668	019010
Calibrator Diluent	101669	019104
Calibrator Concentrate	103217	912801
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	001402	0.588	0.381-0.734	
Control 2	001402	32.5	24.6-36.9	
Panel 1	934301	6.60	1.99-7.85	
Panel 2	934302	24.5	12.1-27.9	
Panel 3	934701	114	72.0-125	

Calibrator Concentrate Value Assignment (pg/mL)	1749
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 21/01/2020

Name

Title

Signature/Date