

Product Number:	101664
Lot Number:	502705
Expiration:	10-Nov-2021
Platform(s):	HD-1/HD-X



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

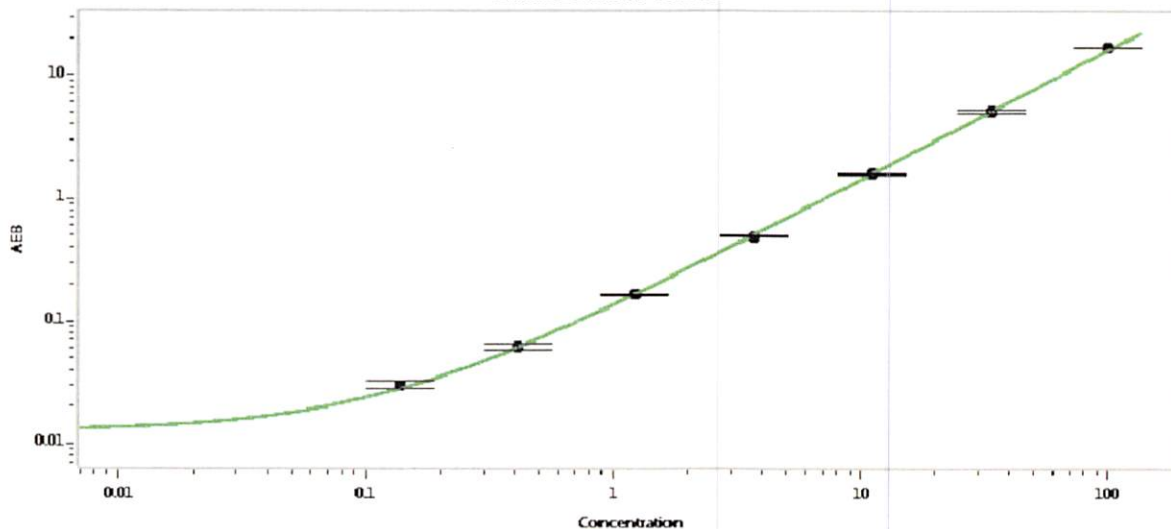
Component	Part Number	Lot Number
Bead Reagent	101666	031516
Detector Reagent	101665	102208
SBG Reagent	101667	032405
Sample Diluent	101668	032415
Calibrator Diluent	101669	032406
Calibrator Concentrate	103217	026801
RGP Reagent	103159	N/A ¹

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	0.905	0.261-1.30	
Control 2	001402	45.2	19.2-62.8	
Panel 1	934301	8.52	3.41-10.1	
Panel 2	934302	31.4	13.2-41.3	
Panel 3	934701	141	58.0-156	

Calibrator Concentrate Value Assignment (pg/mL)	2184
---	------

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

Junpil Park

QC Supervisor

09 Feb 2021

Name

Title

Signature/Date