

Product Number:	103186
Lot Number:	503286
Expiration:	17-May-2022
Platform(s):	HD-1/ HD-X



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

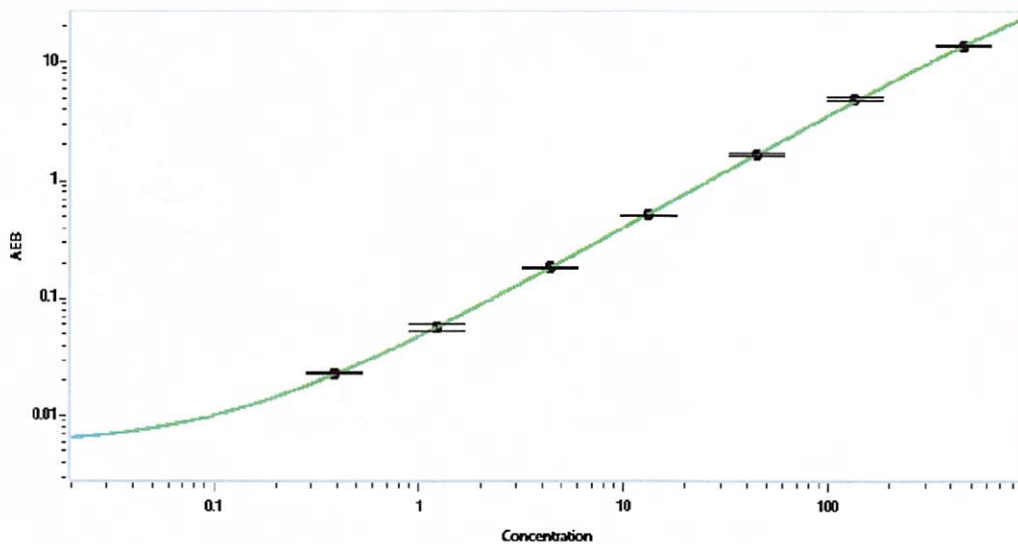
Component	Part Number	Lot Number
Bead Reagent	102246	115708
Detector Reagent	102248	115809
SBG Reagent	102250	115810
Sample Diluent	102252	115914
Calibrators	102255	134704
Control 1	102256	134705
Control 2	102257	134705
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134705	4.01	3.20-4.80
Control 2	134705	271	224-336
Panel 1	134208	6.27	3.95-9.43
Panel 2	134208	3.94	1.86-7.15
Panel 3	134208	69.3	49.9-74.8
Panel 4	135805	366	274-411

* 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	0.391
C	1.22
D	4.35
E	13.2
F	44.8
G	134
H	456

Review/Approval Shivani Goel QC Shivani Goel 27-Jan-2022
 Name Title Signature/Date