

Product Number:	103186
Lot Number:	502637
Expiration:	18-Nov-2021
Platform(s):	HD-1/ HD-X



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

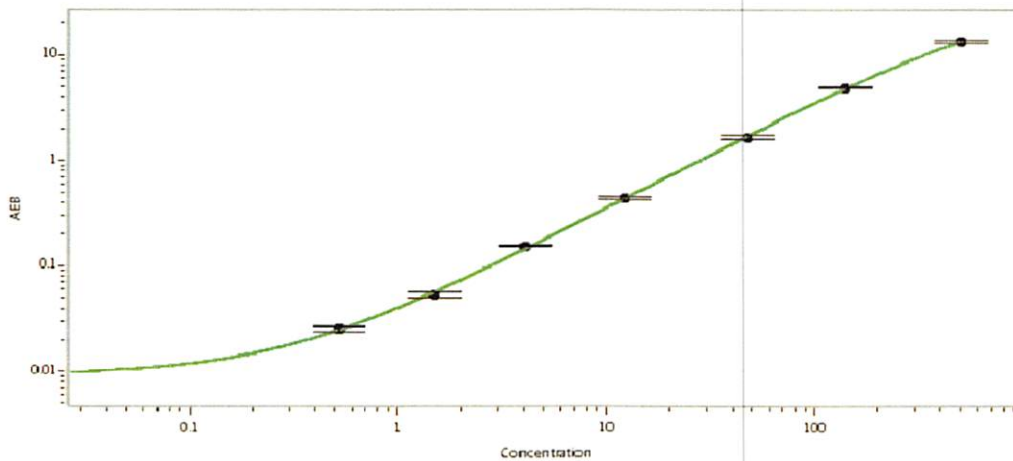
Component	Part Number	Lot Number
Bead Reagent	102246	032812
Detector Reagent	102248	033502
SBG Reagent	102250	033501
Sample Diluent	102252	033505
Calibrators	102255	033513
Control 1	102256	033602
Control 2	102257	033602
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	033602	4.10	3.03-4.59
Control 2	033602	152	125-188
Panel 1	008501	14.7	9.92-18.2
Panel 2	008501	11.2	8.15-13.3
Panel 3	008501	59.7	47.2-84.0
Panel 4	009401	206	144-253

* 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	0.522
C	1.48
D	4.07
E	12.2
F	47.0
G	140
H	496

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

[Signature] 14Dec2020