

Product Number:	103186
Lot Number:	502670
Expiration:	24-Sep-2021
Platform(s):	HD-1/ HD-X



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

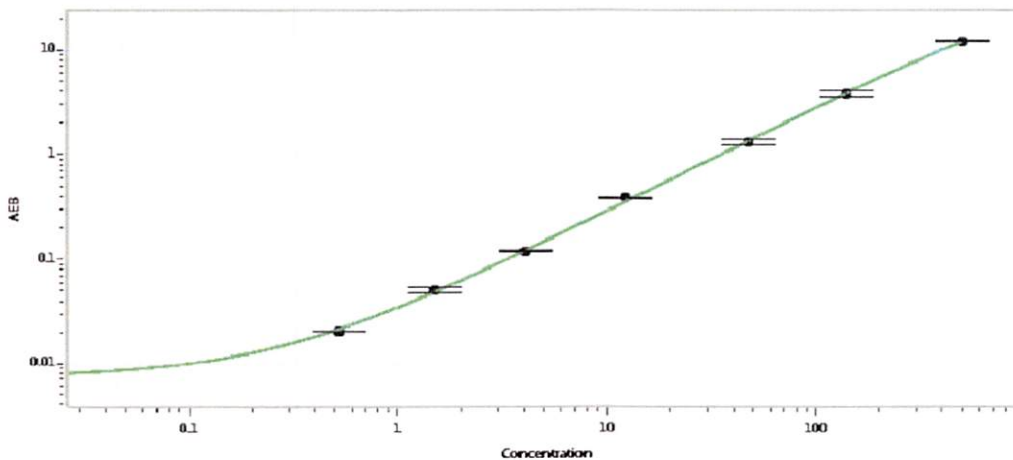
Component	Part Number	Lot Number
Bead Reagent	102246	026804
Detector Reagent	102248	026904
SBG Reagent	102250	026908
Sample Diluent	102252	026905
Calibrators	102255	033513
Control 1	102256	033602
Control 2	102257	033602
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	033602	4.08	2.96-4.77
Control 2	033602	155	125-187
Panel 1	008501	14.5	9.92-18.2
Panel 2	008501	10.6	8.15-13.3
Panel 3	008501	67.0	47.2-84.0
Panel 4	009401	201	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
Name

QC Supervisor
Title

30 Dec 2020
Signature/Date