

Product Number:	103186
Lot Number:	502492
Expiration:	19-Feb-2021
Platform(s):	HD-1/ HD-X



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

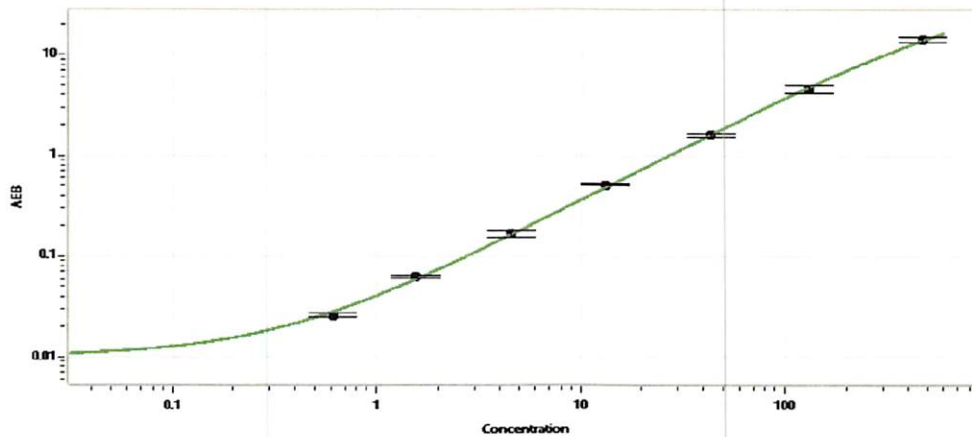
Component	Part Number	Lot Number
Bead Reagent	102246	022302
Detector Reagent	102248	022404
SBG Reagent	102250	022403
Sample Diluent	102252	005002
Calibrators	102255	022401
Control 1	102256	022402
Control 2	102257	022402
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	022402	4.61	3.41-5.21
Control 2	022402	172	132-205
Panel 1	922601	19.9	14.1 - 21.1
Panel 2	922602	6.71	4.94 - 7.41
Panel 3	922603	41.6	28.4 - 42.6
Panel 4	922604	174	129 - 204

\* 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 0.609
- C 1.56
- D 4.56
- E 13.1
- F 43.1
- G 131
- H 468

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 24 Aug 2020

Name

Title

Signature/Date