

Product Number:	103180
Lot Number:	502497
Expiration:	18-Apr-2021
Platform(s):	HD-1/HD-X



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

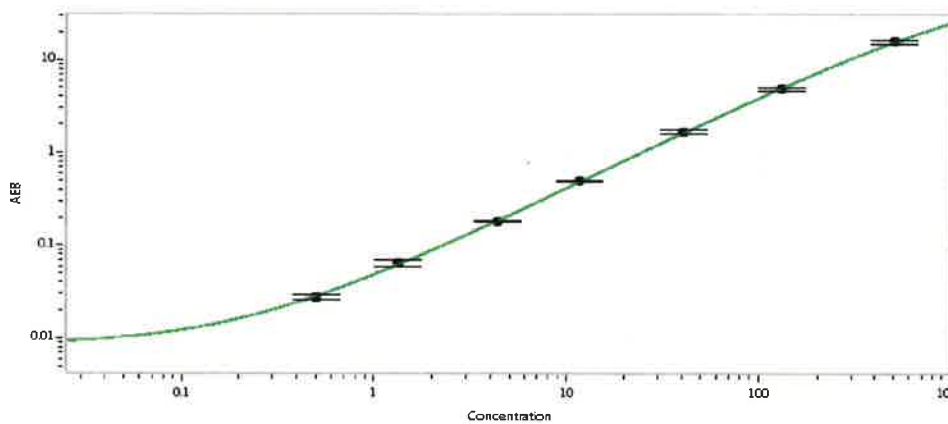
Component	Part Number	Lot Number
Bead Reagent	102246	022302
Detector Reagent	102248	022404
SBG Reagent	102250	022403
Sample Diluent	102252	010903
Calibrators	102255	933601
Control 1	102256	933901
Control 2	102257	933901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	933901	3.51	2.60-4.70
Control 2	933901	158	122-183
Panel 1	922601	18.0	14.1 - 21.1
Panel 2	922602	6.22	4.94 - 7.41
Panel 3	922603	39.2	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 0.501
- C 1.32
- D 4.32
- E 11.8
- F 41.1
- G 131
- H 506

Review/Approval: Shivani Goel Name, QC Manager Title, Shivani Goel 18 Aug 2020 Signature/Date