

Product Number:	103180
Lot Number:	502647
Expiration:	20-Nov-2021
Platform(s):	HD-1/ HD-X



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

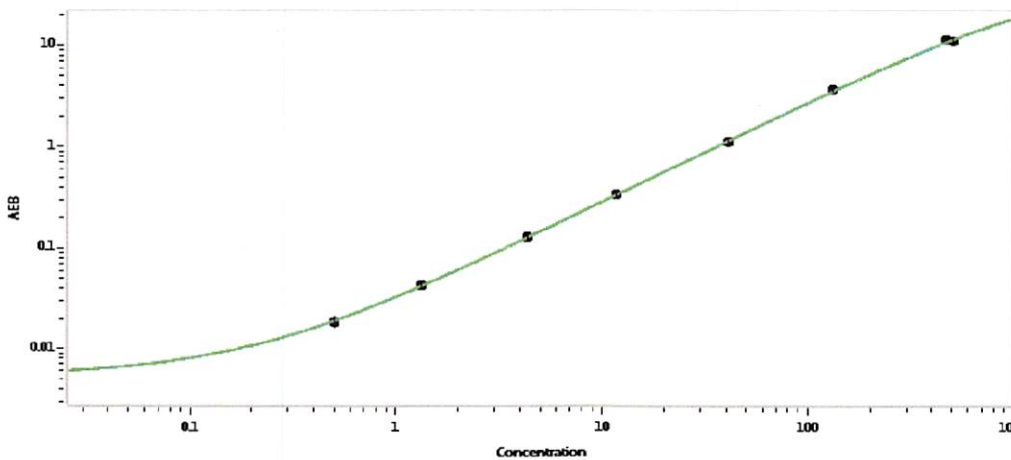
Component	Part Number	Lot Number
Bead Reagent	102246	032812
Detector Reagent	102248	033502
SBG Reagent	102250	033501
Sample Diluent	102252	033505
Calibrators	102255	933601
Control 1	102256	933901
Control 2	102257	933901
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	933901	4.19	2.60 - 4.70
Control 2	933901	178	122 - 183
Panel 1	922601	18.6	14.3 - 21.4
Panel 2	922602	6.62	4.97 - 7.45
Panel 3	922603	38.7	28.8 - 43.2
Panel 4	922604	163	133 - 201

\* 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

Junpil Park  
Name

QC Supervisor  
Title

Signature/Date

*[Handwritten Signature]* 21 Dec 2020