

Product Number:	103186
Lot Number:	503302
Expiration:	21-Jan-2023
Platform(s):	HD-1/ HD-X



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

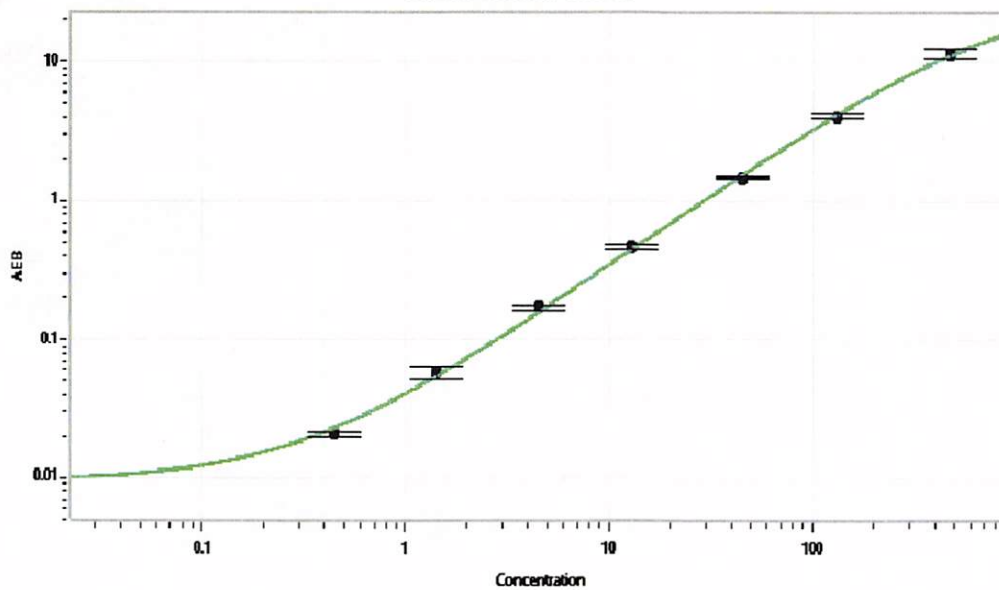
Component	Part Number	Lot Number
Bead Reagent	102246	203806
Detector Reagent	102248	203910
SBG Reagent	102250	203911
Sample Diluent	102252	203912
Calibrators	102255	203104
Control 1	102256	203105
Control 2	102257	203105
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203105	3.53	2.87-4.38
Control 2	203105	196	171-257
Panel 1	134208	7.65	3.95-9.43
Panel 2	134208	4.52	1.86-7.15
Panel 3	134208	62.5	49.9-74.8
Panel 4	135801	303	274-411

\* 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	0.448
C	1.42
D	4.53
E	12.8
F	45.1
G	133
H	472

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 08 Mar 2022

Name

Title

Signature/Date