

Product Number:	103186
Lot Number:	502976
Expiration:	17-May-2022
Platform(s):	HD-1/ HD-X



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

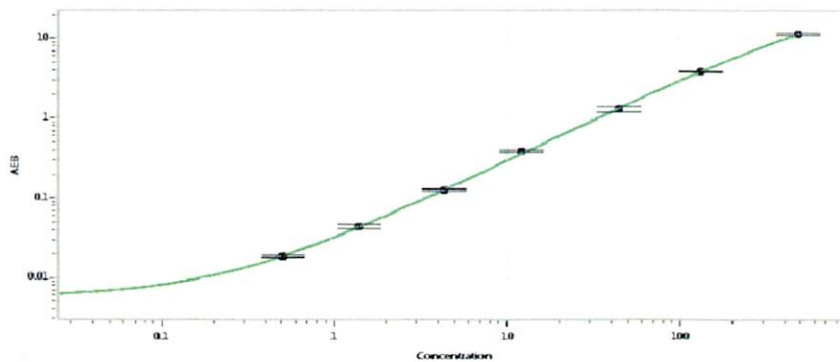
Component	Part Number	Lot Number
Bead Reagent	102246	115708
Detector Reagent	102248	115809
SBG Reagent	102250	115810
Sample Diluent	102252	115914
Calibrators	102255	115906
Control 1	102256	115907
Control 2	102257	115907
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*						* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Previously reported		Corrected **		
		Result (pg/mL)	Mean Range (pg/mL)	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	115907	3.87	3.06-5.03	2.99	2.57-5.23	
Control 2	115907	241	140-314	177	128-312	
Panel 1	008501	15.7	9.26-19.5	11.8	9.26-19.5	
Panel 2	008501	12.1	7.42-14.5	9.15	7.42-14.5	
Panel 3	008501	78.6	46.0-88.2	57.8	46.0-88.2	
Panel 4	009401	275	144-253	201	144-253	

** Values and ranges are updated as of 13Dec2021. Corrected result and calibrator levels should be used for reference.

Calibration Curve



	Calibrator Levels (pg/mL)	
	Previously reported	Corrected**
A	0.000	0.000
B	0.425	0.381
C	1.30	1.02
D	3.97	2.77
E	11.4	8.13
F	38.9	28.7
G	113	89.4
H	364	283

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date 15 Dec 2021