

Product Number:	103186
Lot Number:	502994
Expiration:	28-Jul-2022
Platform(s):	HD-1/ HD-X



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

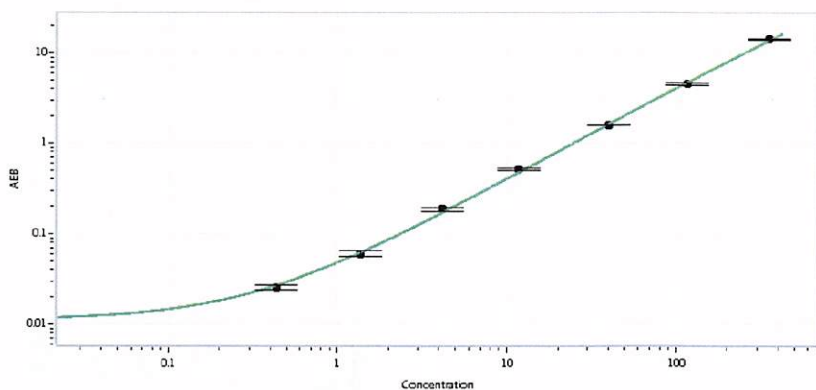
Component	Part Number	Lot Number
Bead Reagent	102246	121015
Detector Reagent	102248	121028
SBG Reagent	102250	121030
Sample Diluent	102252	121031
Calibrators	102255	121411
Control 1	102256	117908
Control 2	102257	117908
RGP Reagent	103159	N/A ¹

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	117908	4.63	3.00-5.12	3.23	2.36-3.55	
Control 2	117908	261	198-297	216	152-228	
Panel 1	008501	17.1	8.59-20.5	12.5	8.59-20.5	
Panel 2	008501	13.4	7.39-15.1	9.69	7.39-15.1	
Panel 3	008501	78.4	43.8-91.8	61.2	43.8-91.8	
Panel 4	009401	249	128-302	206	128-302	

** Values and ranges are updated as of 26Oct2021. 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	0.431	0.269
C	1.36	0.964
D	4.18	3.18
E	11.7	9.49
F	39.7	30.4
G	117	98.5
H	354	340

Review/Approval

Junpil Park
Name

QC Supervisor
Title


Signature/Date
01 Nov 2021