

Product Number:	103186
Lot Number:	502845
Expiration:	30-Mar-2022
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	102246	109016
Detector Reagent	102248	108913
SBG Reagent	102250	108909
Sample Diluent	102252	108914
Calibrators	102255	109806
Control 1	102256	109807
Control 2	102257	109807
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

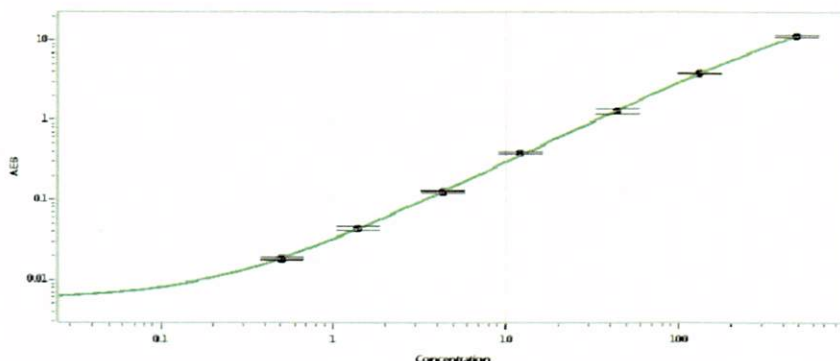


Data below represents results generated on the Simoa HD-1® Analyzer.

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	109807	4.90	3.44-5.87	4.67	3.60 - 5.40	
Control 2	109807	252	174-298	243	183 - 274	
Panel 1	008501	14.1	8.59-20.5	13.6	8.59 - 20.5	
Panel 2	008501	12.0	7.39-15.1	11.5	7.39 - 15.1	
Panel 3	008501	69.3	43.8-91.8	66.5	43.8 - 91.8	
Panel 4	009401	203	144-253	195	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	0.506	0.464
C	1.41	1.38
D	4.33	4.35
E	12.2	11.8
F	44.2	41.5
G	132	126
H	477	507

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date

[Signature] 01 Nov 2021