

Product Number:	102215
Lot Number:	502953
Expiration:	4-Apr-2022
Platform(s):	HD-1/HD-x & SR-X



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

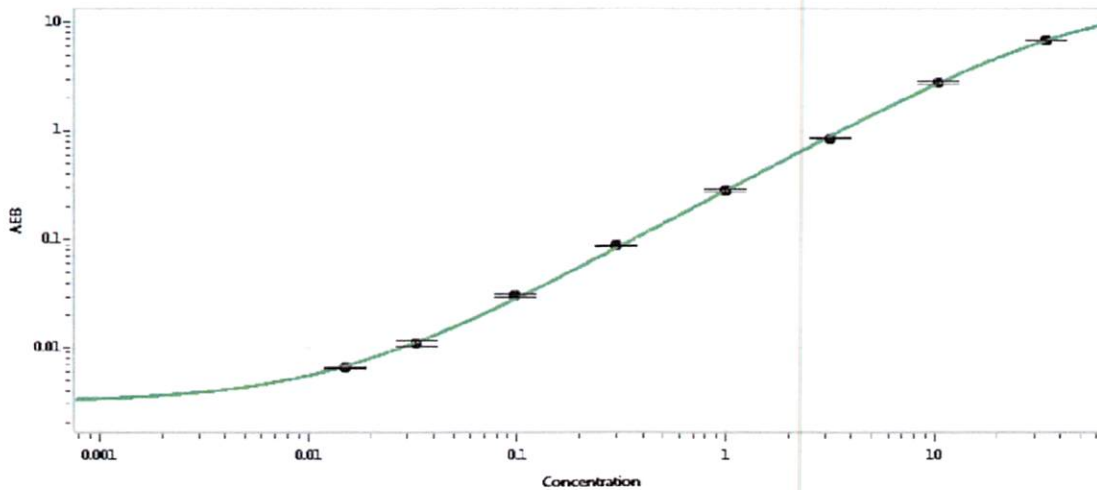
Component	Part Number	Lot Number
Bead Reagent	102211	109401
Detector Reagent	102212	109612
SBG Reagent	102213	109511
Assay Diluent	102214	113910
Calibrators	101310	111111
Control 1	100702	111112
Control 2	100703	111112
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	111112	0.136	0.105-0.157
Control 2	111112	10.6	8.10-13.2
Panel 1	019601	0.196	0.141-0.212
Panel 2	019601	1.87	1.53-2.30
Panel 3	019601	10.7	9.54-14.3

* 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	0.015
C	0.033
D	0.097
E	0.301
F	1.00
G	3.16
H	10.3
I	34.0

Review/Approval Junpil Park QC Supervisor [Signature] 02 Jun 2021
Name Title Signature/Date