

Product Number:	102215
Lot Number:	503088
Expiration:	10-Jul-2022
Platform(s):	HD-1/HD-x & SR-X



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

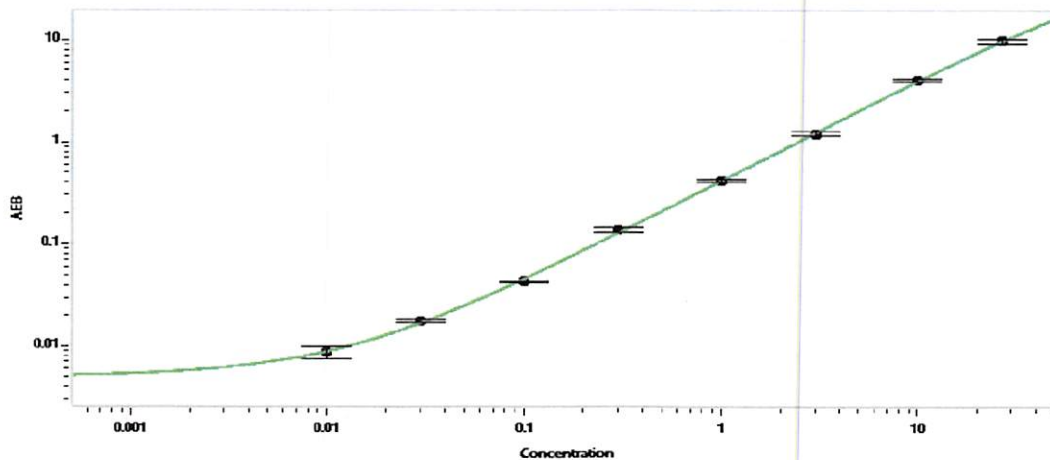
Component	Part Number	Lot Number
Bead Reagent	102211	119101
Detector Reagent	102212	119313
SBG Reagent	102213	119314
Assay Diluent	102214	119905
Calibrators	101310	116506
Control 1	100702	116507
Control 2	100703	116507
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	116507	0.095	0.077 - 0.115
Control 2	116507	9.63	7.25 - 10.9
Panel 1	019601	0.231	0.129 - 0.242
Panel 2	019601	2.26	1.56 - 2.34
Panel 3	019601	12.1	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	0.012
C	0.029
D	0.099
E	0.280
F	0.903
G	2.67
H	9.35
I	28.2

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date

11 Aug 2021