

Product Number:	102215
Lot Number:	502974
Expiration:	17-May-2022
Platform(s):	HD-1/HD-x & SR-X



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

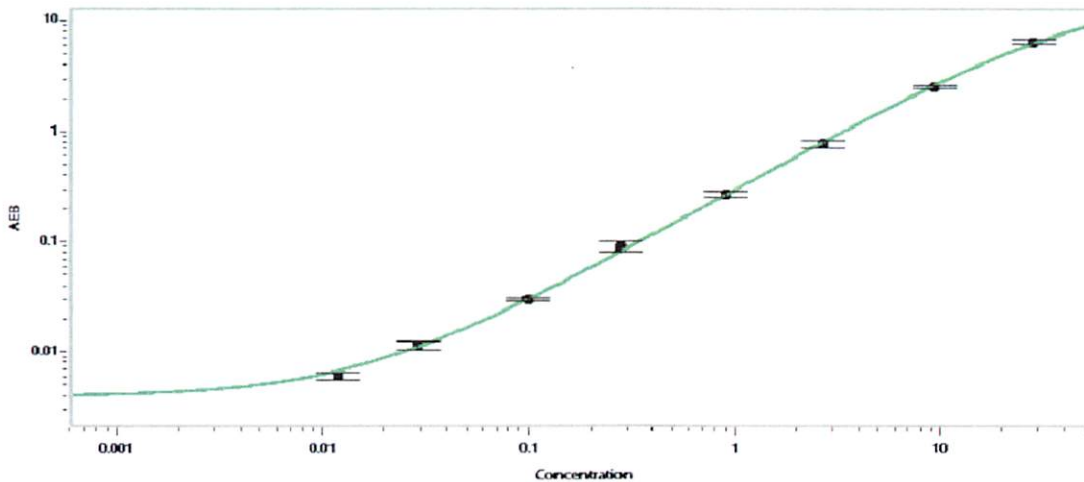
Component	Part Number	Lot Number
Bead Reagent	102211	115601
Detector Reagent	102212	115710
SBG Reagent	102213	115707
Assay Diluent	102214	115711
Calibrators	101310	116506
Control 1	100702	116507
Control 2	100703	116507
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	116507	0.102	0.077-0.115
Control 2	116507	9.24	7.25-10.9
Panel 1	019601	0.206	0.141-0.212
Panel 2	019601	2.02	1.53-2.30
Panel 3	019601	11.8	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

23 Jun 2021