

Product Number:	102215
Lot Number:	502674
Expiration:	29-Dec-2021
Platform(s):	HD-1/HD-x & SR-X



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

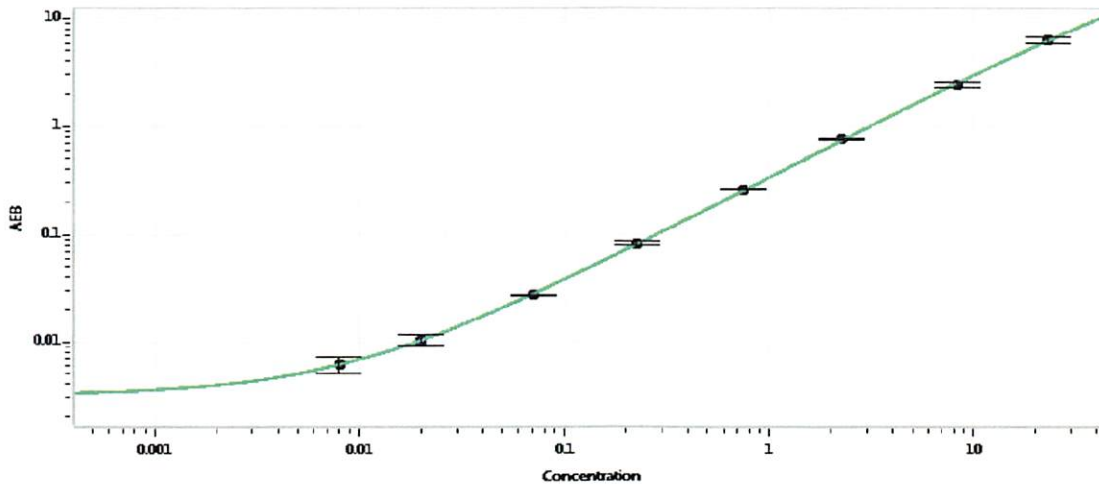
Component	Part Number	Lot Number
Bead Reagent	102211	036409
Detector Reagent	102212	104005
SBG Reagent	102213	104016
Assay Diluent	102214	104912
Calibrators	101310	004401
Control 1	100702	004402
Control 2	100703	004402
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	004402	0.070	0.058-0.087
Control 2	004402	8.14	6.17-9.26
Panel 1	019601	0.180	0.141-0.212
Panel 2	019601	1.99	1.53-2.30
Panel 3	019601	12.4	9.54-14.3

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Review/Approval Junpil Park QC Supervisor [Signature] 03 Mar 2021
Name Title Signature/Date