

Product Number:	102215
Lot Number:	503044
Expiration:	17-Jun-2022
Platform(s):	HD-1/HD-x & SR-X



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

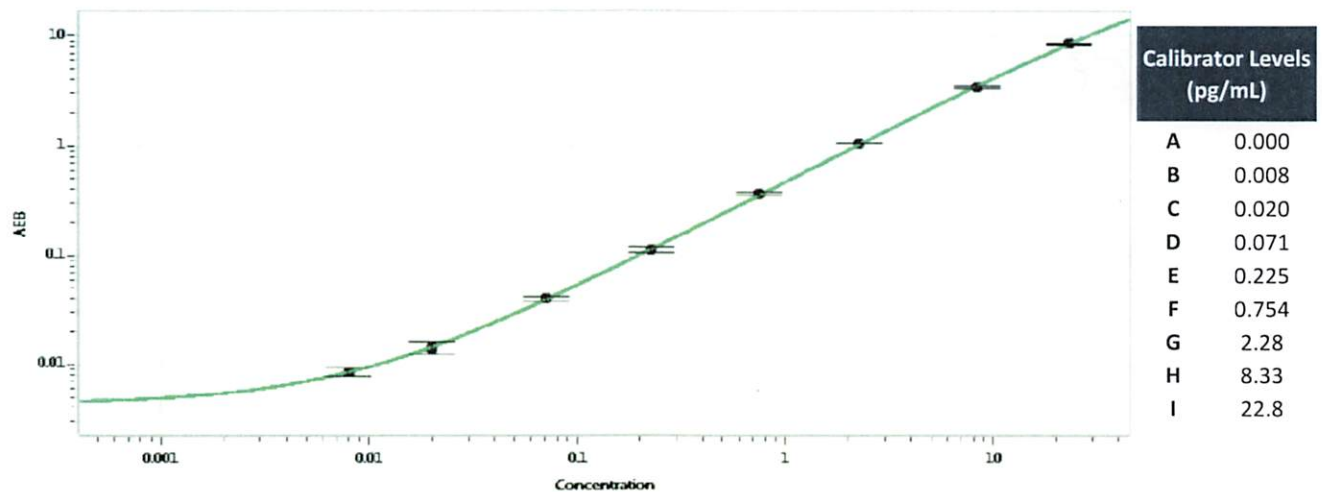
Component	Part Number	Lot Number
Bead Reagent	102211	119101
Detector Reagent	102212	119313
SBG Reagent	102213	119314
Assay Diluent	102214	119905
Calibrators	101310	119903
Control 1	100702	119904
Control 2	100703	119904
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	119904	0.072	0.058-0.126
Control 2	119904	8.61	5.64-9.22
Panel 1	019601	0.175	0.141-0.212
Panel 2	019601	1.97	1.53-2.30
Panel 3	019601	11.9	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 _____
 Name Title Signature/Date
 _____ _____ *[Signature]* 05 Aug 2021