

Product Number:	102215
Lot Number:	502247
Expiration:	30-Oct-2020
Platform(s):	HD-1/HD-x & SR-X



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

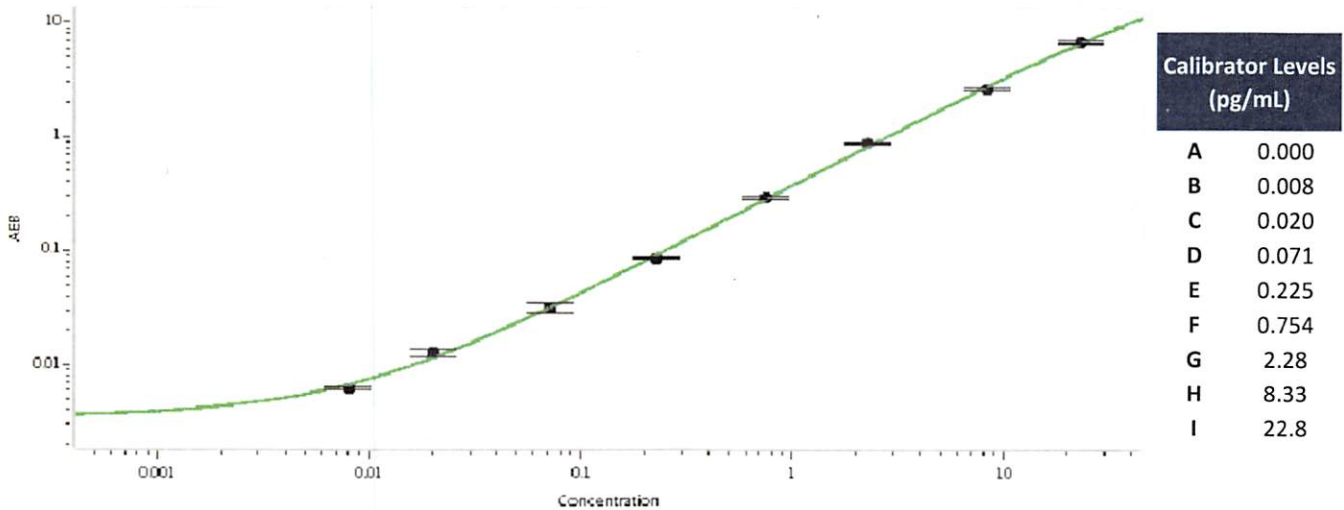
Component	Part Number	Lot Number
Bead Reagent	102211	930810
Detector Reagent	102212	930301
SBG Reagent	102213	930302
Sample Diluent	102214	930303
Calibrators	101310	004401
Controls	100702-100703	004402
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	004402	0.072	0.057-0.086
Control 2	004402	7.48	6.29-9.44
Panel 1	MF-092217-01	0.282	0.177-0.353
Panel 2	MF-092217-02	3.01	1.99-3.51
Panel 3	MF-101117-01	19.0	12.8-20.8

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

Calibration Curve



Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stadler 27 Feb 2020
 Name Title Signature/Date