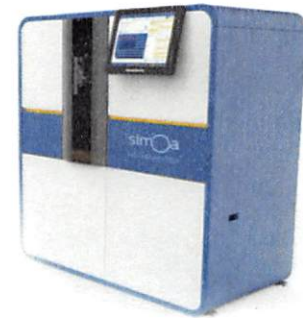


Product Number:	103186
Lot Number:	501912
Expiration:	16-Feb-2020
Platform(s):	HD-1/HD-x & SR-X



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

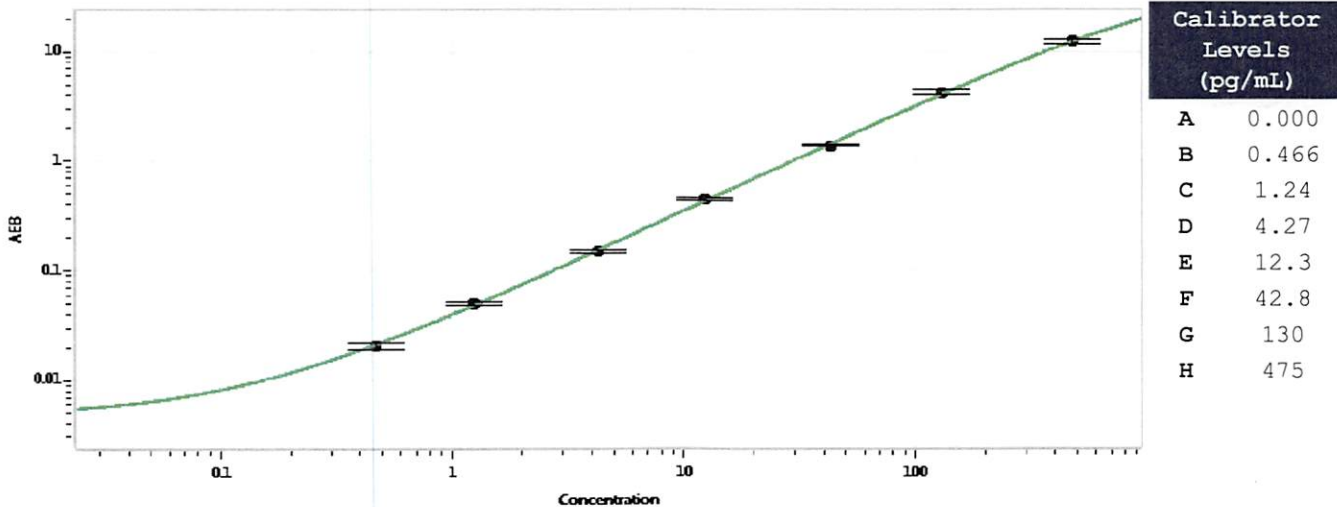
Component	Part Number	Lot Number
Bead Reagent	102246	916103
Detector Reagent	102248	913005
SBG Reagent	102250	913001
Sample Diluent	102252	904701
Calibrators	102255	918901
Controls	102256 - 102257	918902
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918902	3.49	2.28-4.03
Control 2	918902	135	102-152
Panel 1 plasma	823301	5.04	4.08-6.68
Panel 1 serum	823302	6.57	4.90-7.61
Panel 2	823601	80.8	59.3-101
Panel 3	823602	298	236-402

* 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 17 Jul 2019
 Name Title Signature/Date