

Product Number:	103186
Lot Number:	501867
Expiration:	18-Apr-2020
Platform(s):	HD-1, HD-x & SR-X



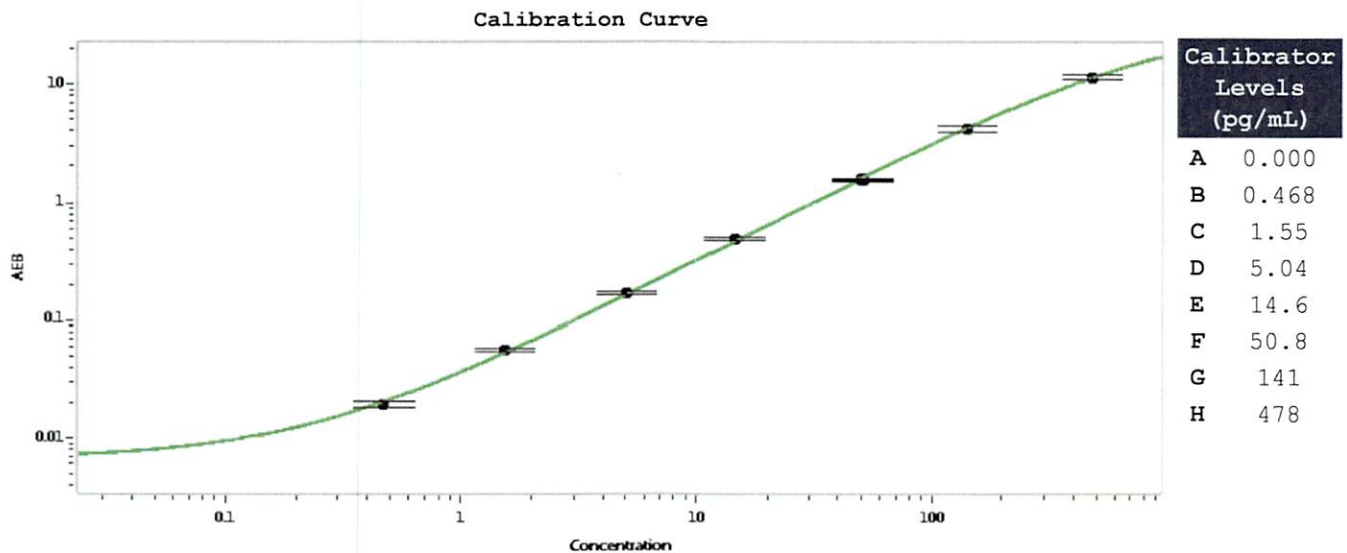
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	916103
Detector Reagent	102248	913005
SBG Reagent	102250	913001
Sample Diluent	102252	910803
Calibrators	102255	916301
Controls	102256 - 102257	916401
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	916401	4.44	3.30-5.02
Control 2	916401	169	133-199
Panel 1a	823301	5.23	4.08-6.68
Panel 1b	823302	6.57	4.90-7.61
Panel 2	823601	80.6	59.3-101
Panel 3	823602	326	236-402

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



Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 24 JUN 2019
	Name	Title	Signature/Date