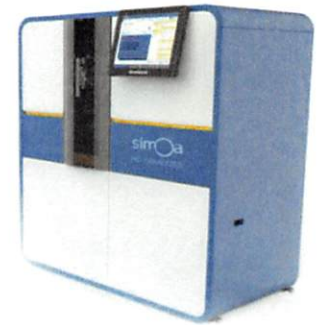


Product Number:	102215
Lot Number:	501949
Expiration:	17-Jan-2020
Platform(s):	HD-1/ HD-X & SR-X



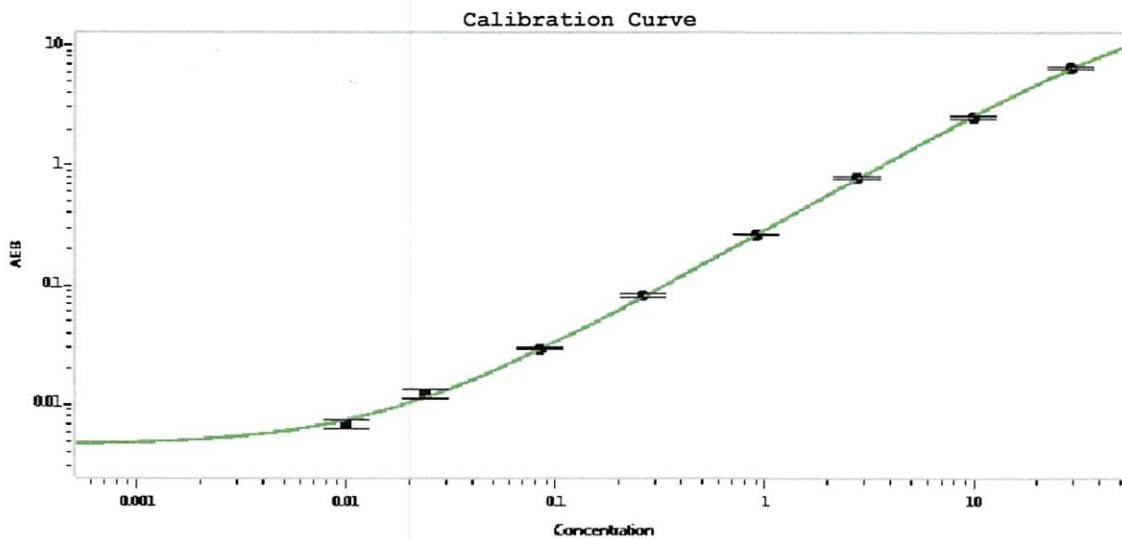
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102211	917007
Detector Reagent	102212	901704
SBG Reagent	102213	916815
Sample Diluent	102214	916814
Calibrators	101310	917801/919005
Controls	100702-100703	917802
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	917802	0.100	0.076-0.114
Control 2	917802	10.8	8.40-12.6
Panel 1	MF-092217-01	0.282	0.177-0.353
Panel 2	MF-092217-02	2.69	1.99-3.51
Panel 3	MF-101117-01	16.9	12.8-20.8

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



Calibrator Levels (pg/mL)

A	0.000
B	0.010
C	0.024
D	0.085
E	0.261
F	0.908
G	2.75
H	9.85
I	28.7

Review/Approval: Patrizia Stadler Assoc. Director, QC *P Stadler* 30 Jul 2019
Name Title Signature/Date