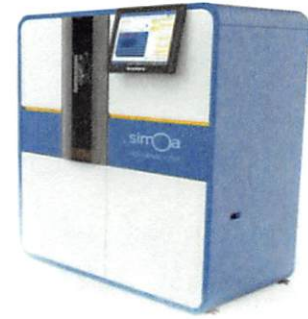


Product Number:	103186
Lot Number:	501953
Expiration:	1-Jul-2020
Platform(s):	HD-1/HD-x & SR-X



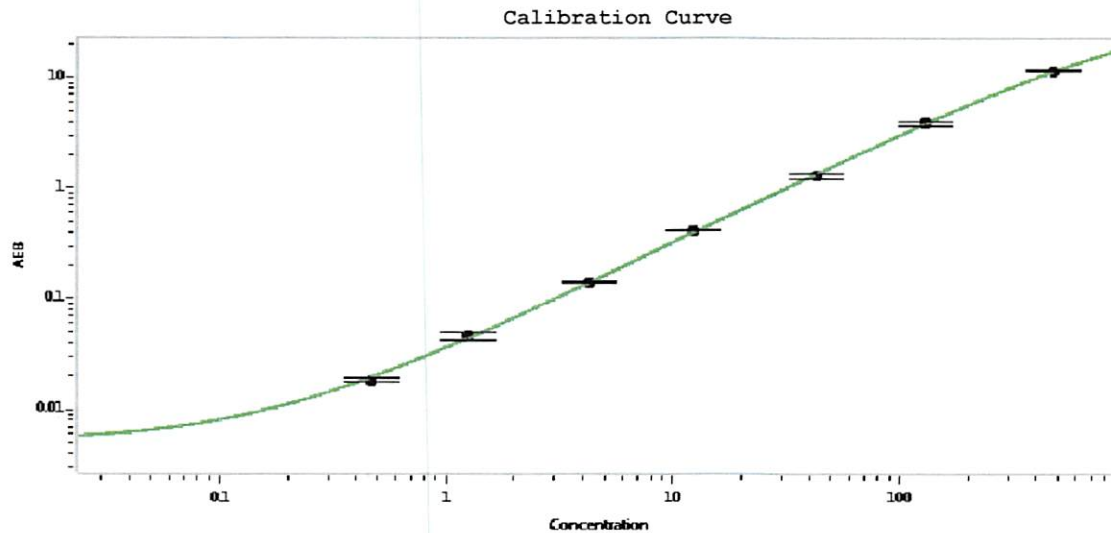
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	920506
Detector Reagent	102248	918213
SBG Reagent	102250	920705
Sample Diluent	102252	919214
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	3.41	2.30-4.01
Control 2	918902	124	102-154
Panel 1a	823301	4.71	4.08-6.68
Panel 1b	823302	6.19	4.90-7.61
Panel 2	823601	79.7	59.3-101
Panel 3	823602	318	236-402

* 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	0.466
C	1.24
D	4.27
E	12.3
F	42.8
G	130
H	475

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	30 Jul 2019
	Name	Title	Signature/Date	