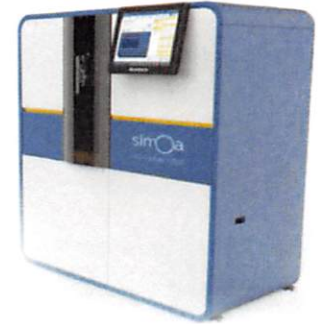


Product Number:	103186
Lot Number:	501683
Expiration:	15-Feb-2020
Platform(s):	HD-1 & SR-X



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

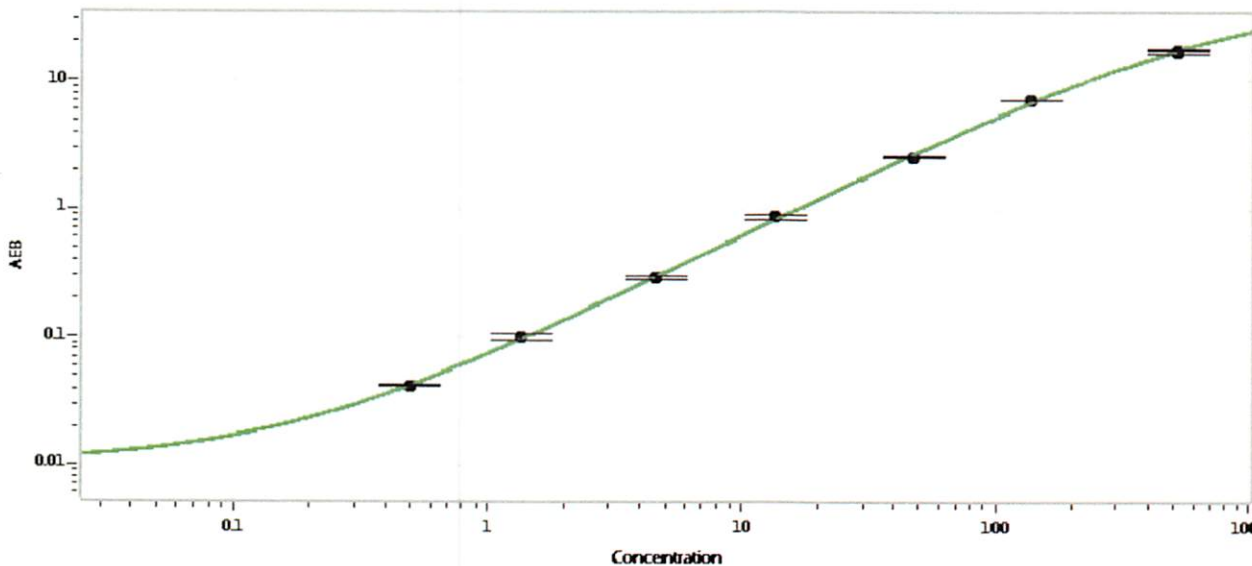
Component	Part Number	Lot Number
Bead Reagent	102246	905611
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Reference Calibrators	102255	905701
Controls	102256 - 102257	905702
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	903502	3.86	3.25-4.87
Control 2	903502	170	134-200
Panel 1-A	823301	5.49	4.08-6.68
Panel 1-B	823302	5.75	4.90-7.61
Panel 2	823601	83.2	59.3-101
Panel 3	823602	330	236-402

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.499
C	1.36
D	4.63
E	13.6
F	47.2
G	136
H	520

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	04 Mar 2019
	Name	Title	Signature/Date	