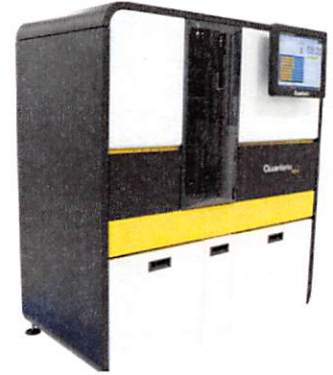


Product Number:	103186
Lot Number:	502031
Expiration:	5-Dec-2019
Platform(s):	HD-1/HD-x & SR-X



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

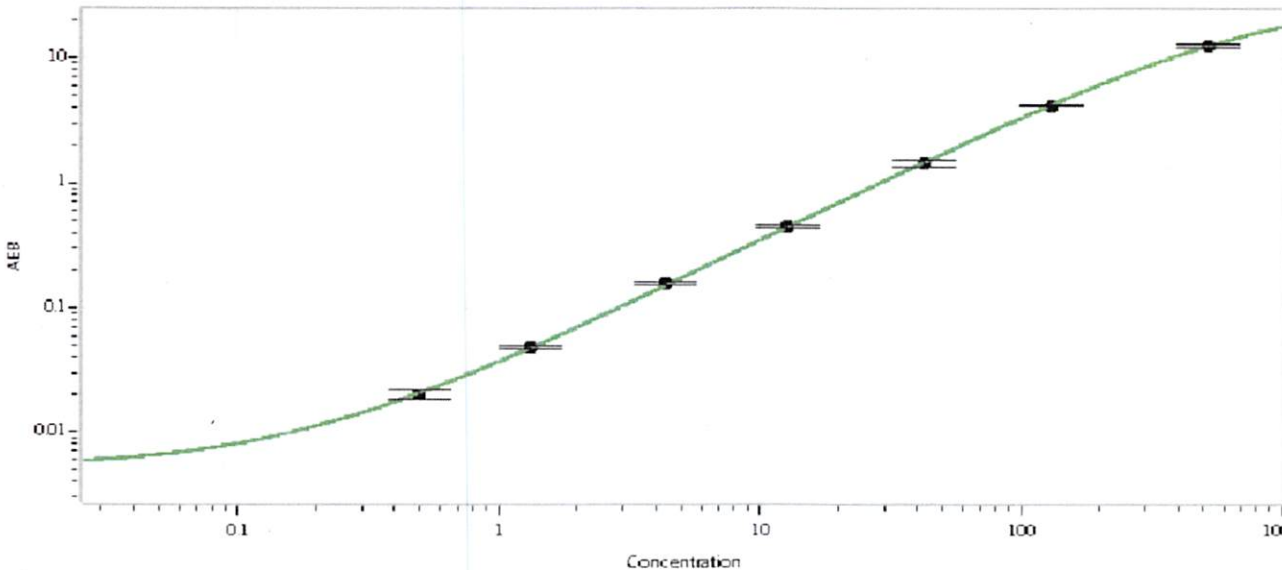
Component	Part Number	Lot Number
Bead Reagent	102246	916103
Detector Reagent	102248	834705
SBG Reagent	102250	833911
Sample Diluent	102252	834507
Calibrators	102255	924701
Controls	102256 - 102257	924803
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	924803	4.21	2.50-4.49
Control 2	924803	135	116-173
Panel 1a	823301	4.58	4.08-6.68
Panel 1b	823302	5.82	4.90-7.61
Panel 2	823601	74.2	59.3-101
Panel 3	823602	283	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	0.503
C	1.33
D	4.32
E	12.8
F	42.3
G	130
H	521

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p. Stadler</i>
	Name	Title	Signature/Date
			18 Sep 2019