

Product Number:	103186
Lot Number:	501733
Expiration:	16-Feb-2020
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	906503
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Reference Calibrators	102255	907801
Controls	102256 - 102257	907802
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

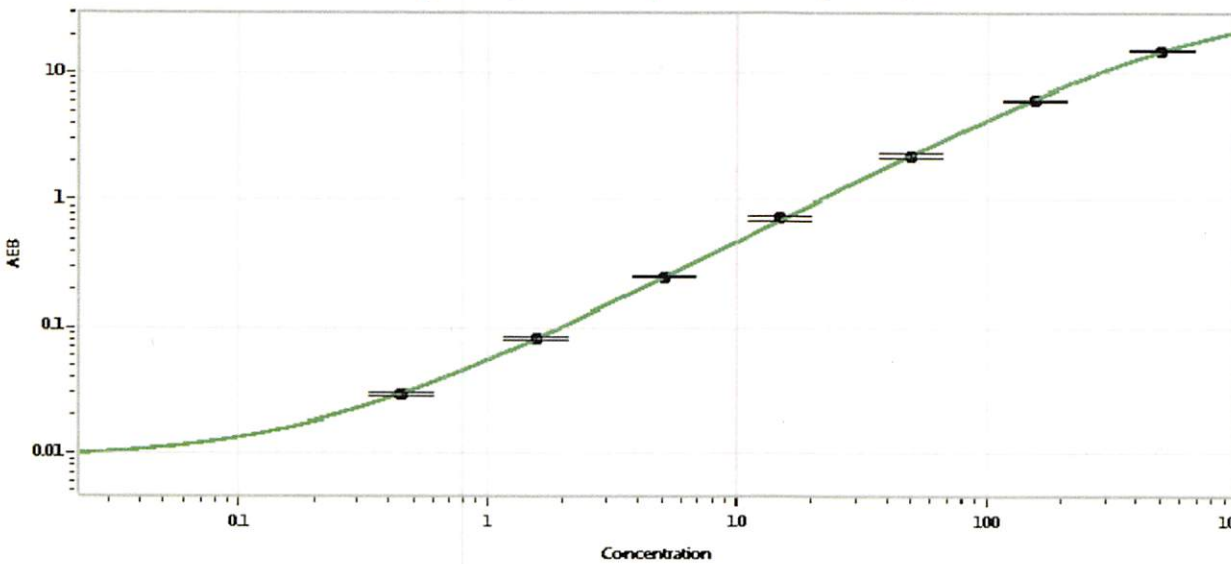


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	907802	4.39	3.25-4.88
Control 2	907802	162	128-192
Panel 1-A	823301	4.98	4.08-6.68
Panel 1-B	823302	5.78	4.90-7.61
Panel 2	823601	80.1	59.3-101
Panel 3	823602	286	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	0.450
C	1.56
D	5.12
E	14.8
F	49.6
G	156
H	504

Review/Approval

Patrizia Stadler QC Assoc. Director
Name Title

P. Stadler 26 MAR 2019
Signature/Date