

Product Number:	103186
Lot Number:	501769
Expiration:	3-Apr-2020
Platform(s):	HD-1 & SR-X



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

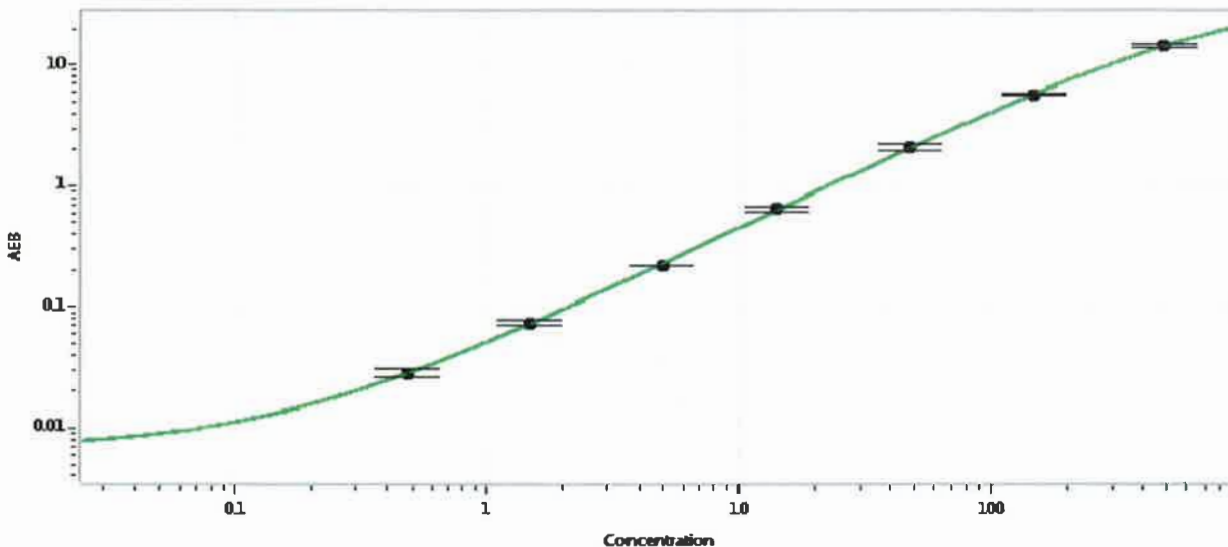
Component	Part Number	Lot Number
Bead Reagent	102246	909315
Detector Reagent	102248	910008
SBG Reagent	102250	909314
Sample Diluent	102252	909302
Reference Calibrators	102255	909905
Controls	102256 - 102257	909906
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909906	3.72	3.22-4.97
Control 2	909906	155	137-205
Panel 1-A	823301	5.03	4.08-6.68
Panel 1-B	823302	5.61	4.90-7.61
Panel 2	823601	75.0	59.3-101
Panel 3	823602	296	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	0.480
C	1.47
D	4.93
E	14.1
F	47.8
G	149
H	482

Review/Approval

Patrizia Stadler QC Assoc. Director
Name Title

P. Stadler 18 Apr 2019
Signature/Date