

Product Number:	103186
Lot Number:	501705
Expiration:	16-Feb-2020
Platform(s):	HD-1 & SR-X



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

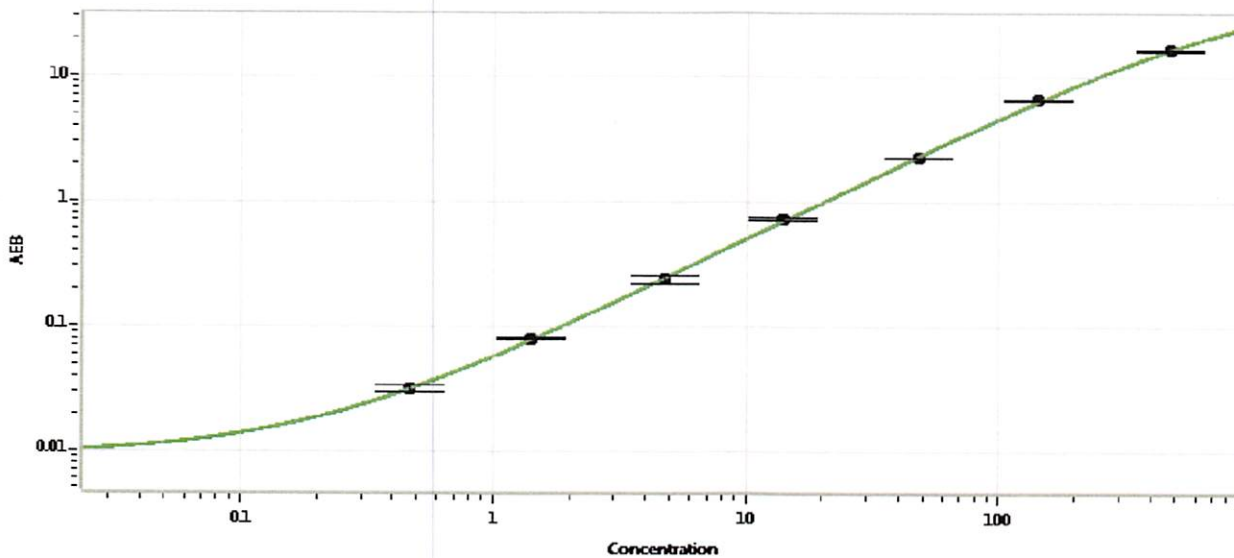
Component	Part Number	Lot Number
Bead Reagent	102246	905611
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Reference Calibrators	102255	906401
Controls	102256 - 102257	906402
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	906402	4.37	3.06-4.94
Control 2	906402	161	129-193
Panel 1-A	823301	5.44	4.08-6.68
Panel 1-B	823302	6.64	4.90-7.61
Panel 2	823601	88.9	59.3-101
Panel 3	823602	313	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	0.466
C	1.42
D	4.79
E	14.0
F	47.6
G	142
H	472

Review/Approval: Patrizia Stadler QC Assoc. Director P. Stadler 11/16/2019  
Name Title Signature/Date