

Product Number:	101622
Lot Number:	501580
Expiration:	19-Sep-2019
Platform(s):	HD-1 & SR-X



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

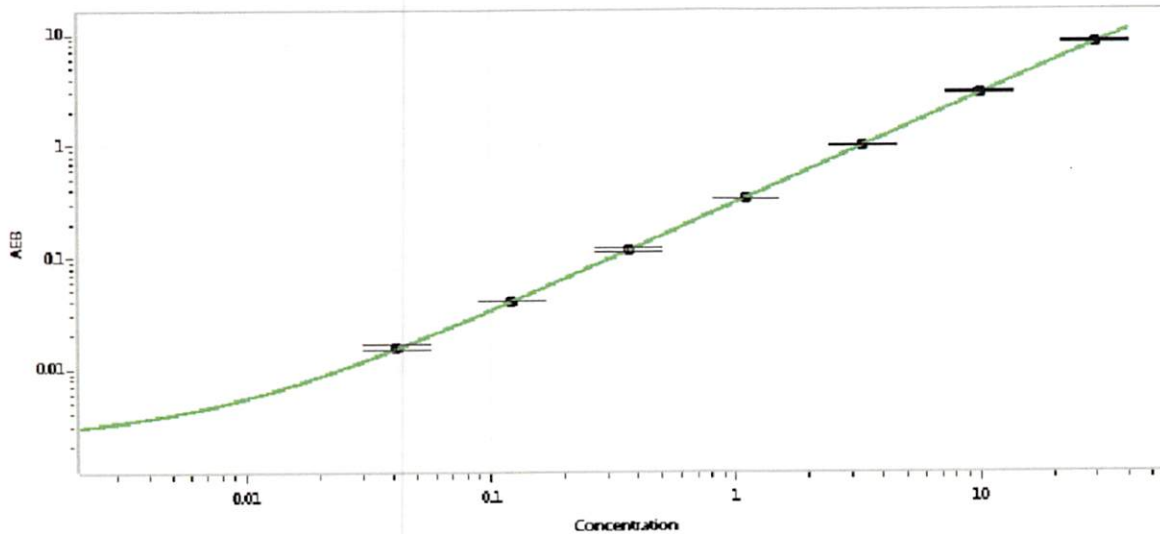
Component	Part Number	Lot Number
Bead Reagent	101624	900918
Detector Reagent	101623	900803
SBG Reagent	101625	900705
Sample Diluent	103203	826225
Calibrator Diluent	101627	834711
Calibrator Concentrate	100652	72350
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	52173	1.84	1.27-1.91	
Control 2	52173	29.6	19.1-31.4	
Panel 1	MF-050918-01	58.7	41.3-66.4	
Panel 2	MF-050918-02	7.88	5.08-8.55	
Panel 3	MF-050918-03	1.02	0.635-1.22	

Calibrator Concentrate Value Assignment (pg/mL)	797
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i> 18 JAN 2019
	Name	Title	Signature/Date