

Product Number:	101605
Lot Number:	501838
Expiration:	27-Feb-2020
Platform(s):	HD-1



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

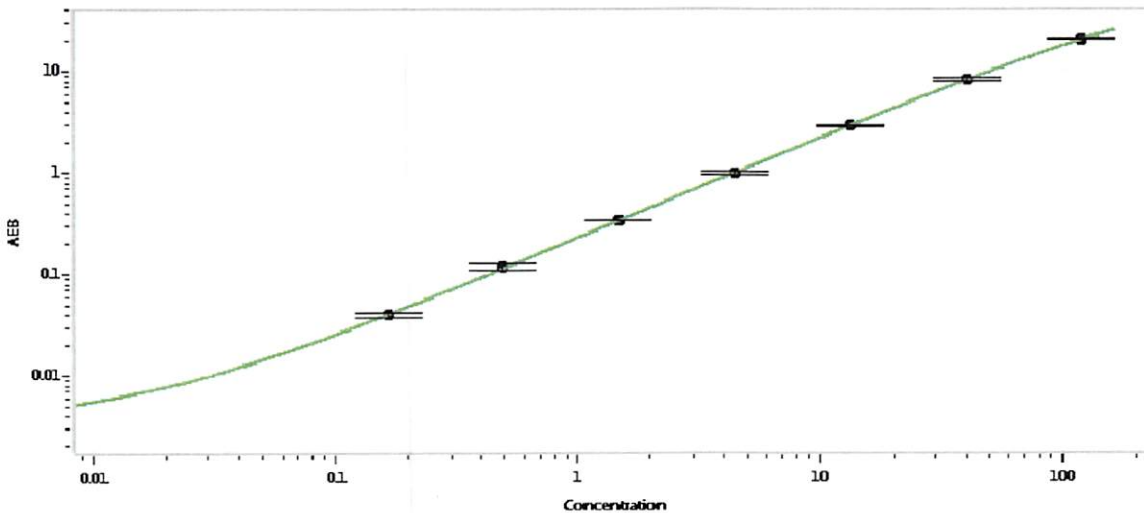
Component	Part Number	Lot Number
Bead Reagent	101607	905810
Detector Reagent	101606	906403
SBG Reagent	101608	906304
Sample Diluent	103200	906309
Calibrator Diluent	101610	906014
Calibrator Concentrate	100895	912609
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	911406	2.37	1.18-2.76	
Control 2	911406	145	81.3-190	
Panel 1	829001	2.76	1.85-4.32	
Panel 2	829002	2.48	2.17-5.07	
Panel 3	829003	9.20	6.71-15.7	
Panel 4	829004	48.6	40.7-61.1	

Calibrator Concentrate Value Assignment (pg/mL)	6,163
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval: Patrizia Stadler Associate Director, QC  
 Name: Patrizia Stadler Title: Associate Director, QC Signature/Date: P. Stadler 2017y 2019