

Product Number:	101672
Lot Number:	501794
Expiration:	29-Jan-2020
Platform(s):	HD-1



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

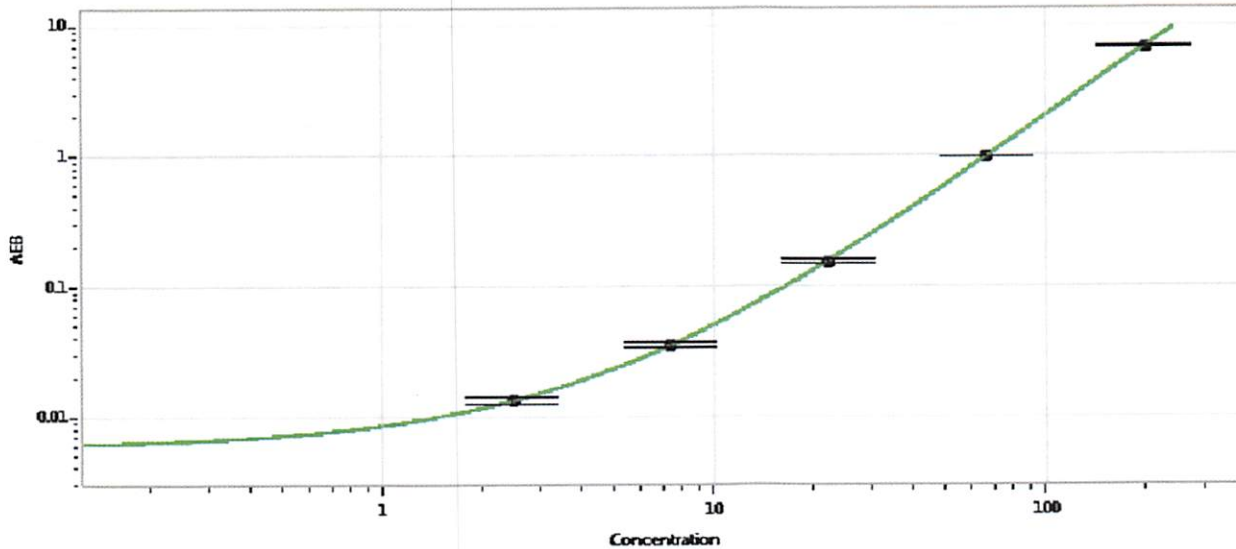
Component	Part Number	Lot Number
Bead Reagent	101674	906011
Detector Reagent	101673	905801
SBG Reagent	101675	905803
Sample Diluent	101676	902929
Calibrator Diluent	101677	904415
Calibrator Concentrate	101734	911207
RGP Reagent	103159	N/A ¹

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	MF-021816-02	22.4	17.1-28.3	
Control 2	MF-021816-02	300	263-434	
Panel 1	MF-040518-01	150	138-207	
Panel 2	MF-040518-02	41.2	33.8-60.0	

Calibrator Concentrate Value Assignment (pg/mL)	5847
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200

Review/Approval	Patrizia Stadler	QC Assoc Director	<i>p Stadler</i>	25 Apr 2019
	Name	Title	Signature/Date	