

Product Number:	101154
Lot Number:	502237
Expiration:	12-Apr-2020
Platform(s):	HD-1/HD-x



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

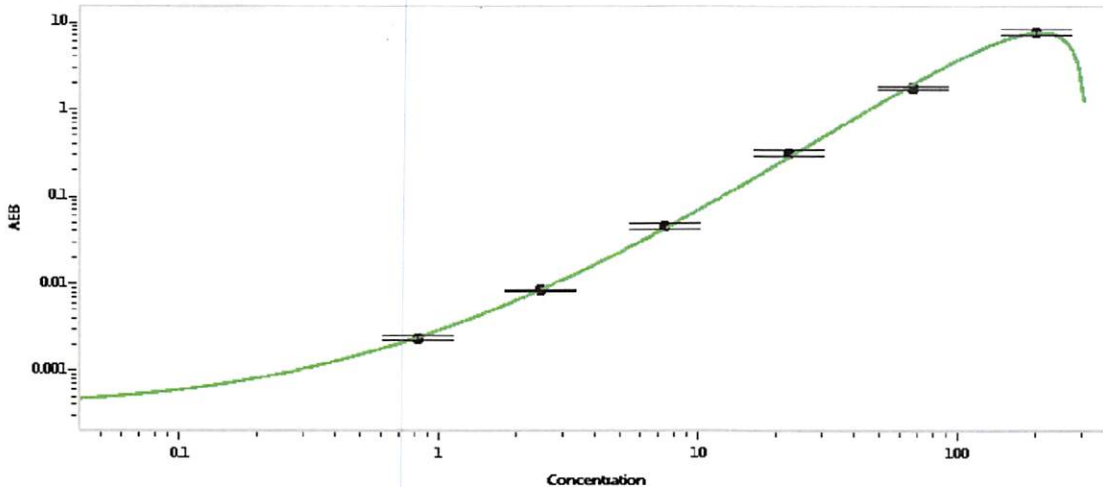
Component	Part Number	Lot Number
Bead Reagent	101141	004510
Detector Reagent	101143	004108
SBG Reagent	101153	004109
Sample Diluent	103320	933816
Calibrator Diluent	101147	934513
Calibrator Concentrate	101145	004202
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918402	43.8	31.9-47.9
Control 2	918402	402	248-453
Panel 1	909901	91.7	68.5-103
Panel 2	909902	77.8	61.9-93.4
Panel 3	909903	295	166-336
Panel 4	909904	288	166-314

* 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	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	26 Feb 2020
	Name	Title	Signature/Date	