

Product Number:	100988
Lot Number:	502145
Expiration:	4-Oct-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	100974	932915
Detector Reagent	100976	932913
SBG Reagent	100987	932701
Sample Diluent	100985	927724
Calibrator Diluent	100980	932620
Calibrator Concentrate	100978	934402
RGP Reagent	103159	N/A ¹

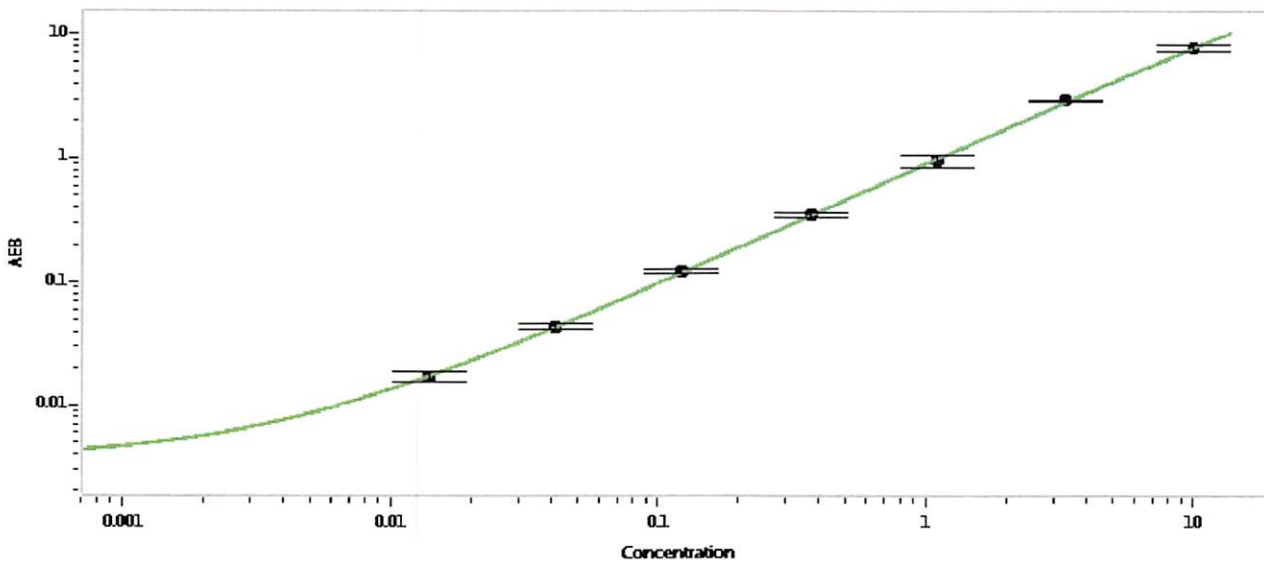
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-101816-02	0.110	0.068-0.161
Control 2	MF-101816-02	9.15	7.32-11.2
Panel 1	MF-101716-04	0.242	0.167-0.327
Panel 2	MF-101716-05	0.062	0.021-0.087
Panel 3	MF-101816-03	18.6	15.6-23.9
Panel 4	MF-050118-01	0.221	0.150-0.293

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	225
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	03 Jan 2020
	Name	Title	Signature/Date	