

Product Number:	100988
Lot Number:	501896
Expiration:	21-Dec-2019
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	910009
Detector Reagent	100976	908007
SBG Reagent	100987	908006
Sample Diluent	100985	907306
Calibrators	100980	835508
Controls	100978	828407
RGP Reagent	103159	N/A ¹

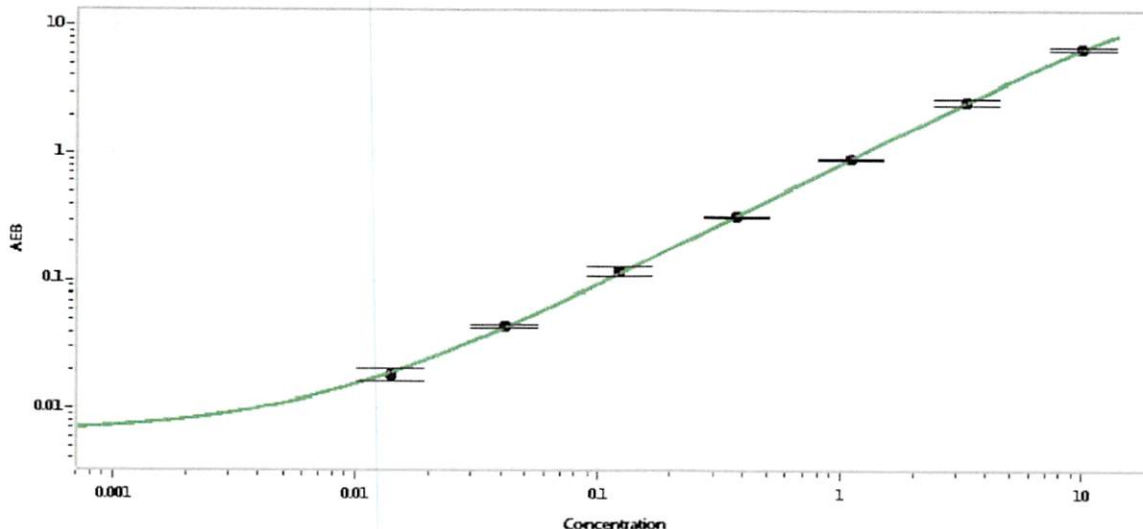
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-101816-02	0.112	0.066-0.163
Control 2	MF-101816-02	9.40	7.28-11.0
Panel 1	MF-101716-04	0.259	0.167-0.317
Panel 2	MF-101716-05	0.053	0.022-0.081
Panel 3	MF-101816-03	20.3	15.6-23.4
Panel 4	MF-050118-01	0.214	0.148-0.287

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

Calibrator Concentrate Value Assignment (pg/mL)	243
---	-----

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>	17 Jul 2019
	Name	Title	Signature/Date	