

Product Number:	101672
Lot Number:	502055
Expiration:	30-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	101674	926106
Detector Reagent	101673	926005
SBG Reagent	101675	926002
Sample Diluent	101676	912026
Calibrator Diluent	101677	924906
Calibrator Concentrate	101734	927002
RGP Reagent	103159	N/A <sup>1</sup>

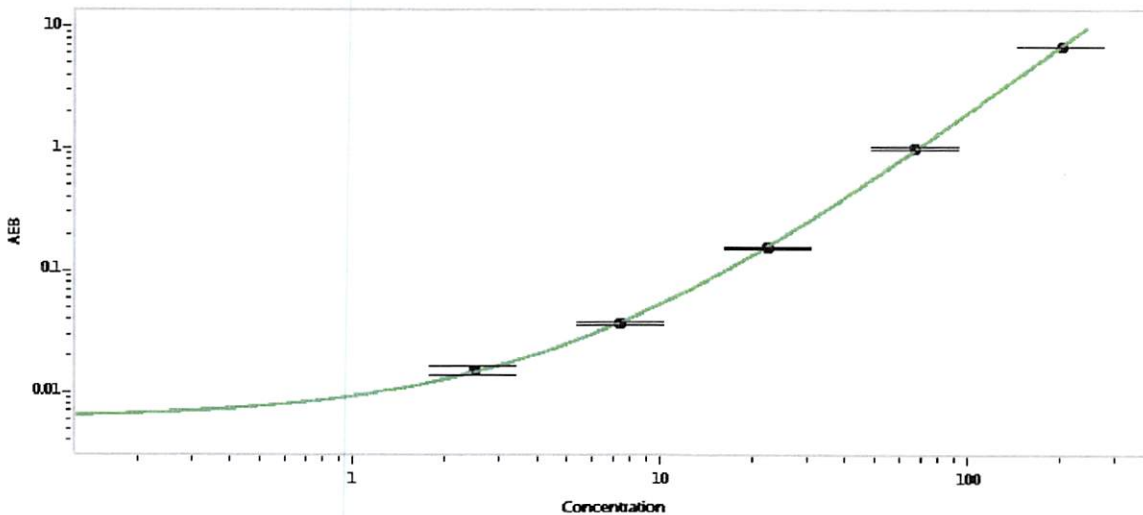
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918301	15.0	11.5-18.9
Control 2	919901	414	325-487
Panel 1	MF-040518-01	179	138-207
Panel 2	MF-040518-02	54.2	33.8-60.0

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

Calibrator Concentrate Value Assignment (pg/mL)	4991
---	------

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

*P. Stadler*

24 Oct 2019

Name

Title

Signature/Date