

Product Number:	101672
Lot Number:	502130
Expiration:	17-Sep-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	927720
Calibrator Diluent	101677	930818
Calibrator Concentrate	101734	927002
RGP Reagent	103159	N/A ¹

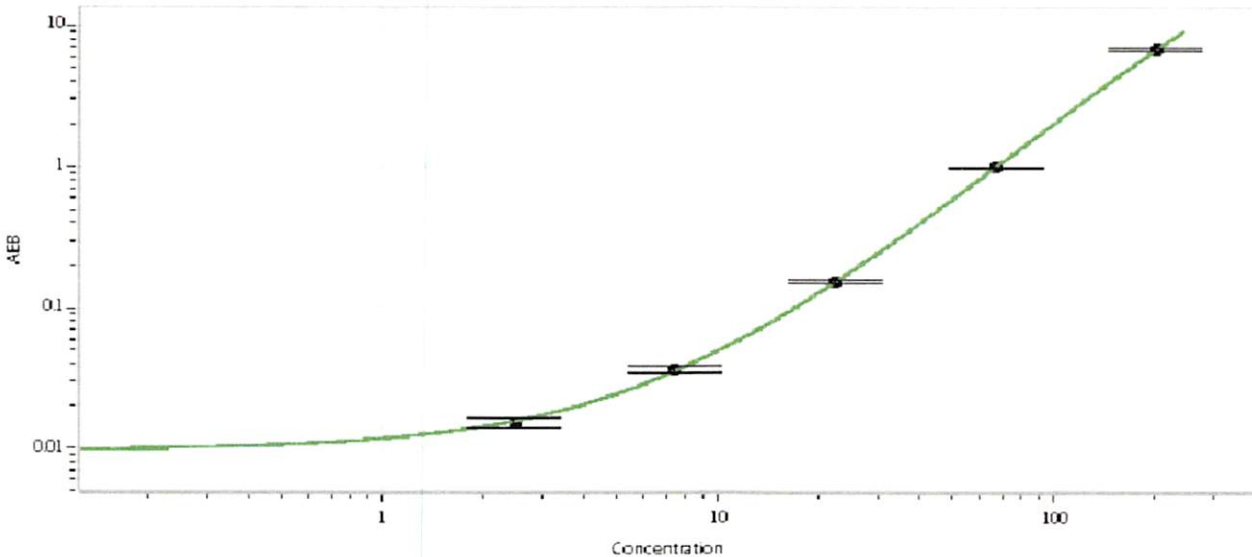
¹ RGP is not Kit Lot Specific

	Release Materials*		
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918301	18.0	10.8-20.7
Control 2	919901	437	330-495
Panel 1	MF-040518-01	201	144-215
Panel 2	MF-040518-02	62.3	30.4-69.0

* 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
Name

QC Assoc Director
Title

P. Stadler 09 Dec 2019
Signature/Date