

Product Number:	101672
Lot Number:	502452
Expiration:	21-Feb-2021
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	017705
Detector Reagent	101673	014205
SBG Reagent	101675	014208
Sample Diluent	101676	005232
Calibrator Diluent	101677	012307
Calibrator Concentrate	101734	911207
RGP Reagent	103159	N/A ¹

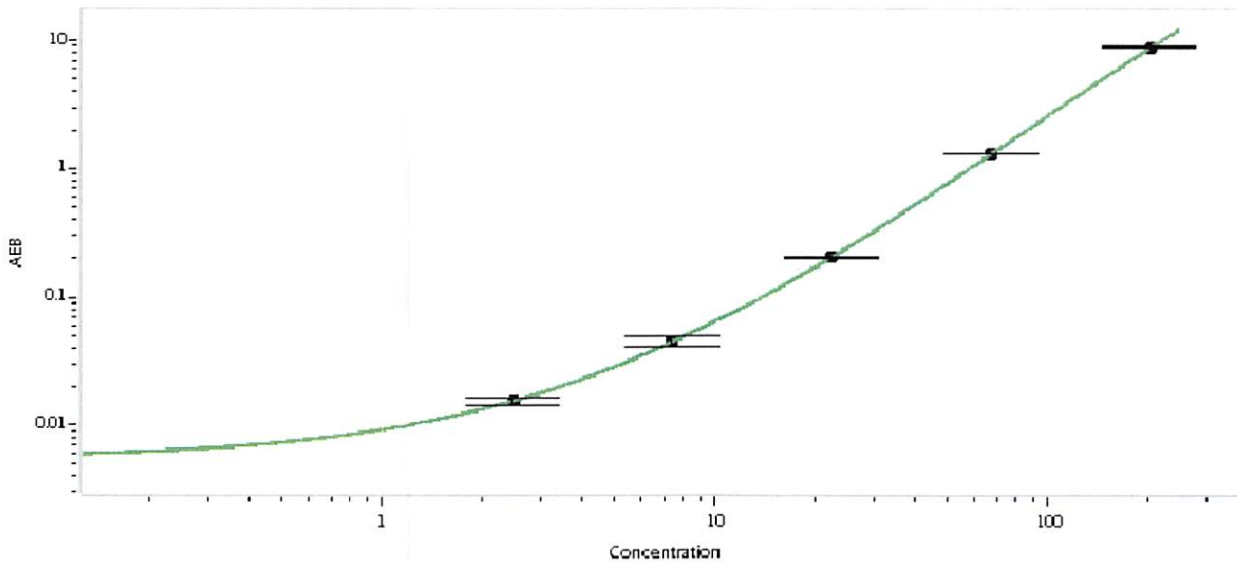
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918301	15.8	10.8-20.7
Control 2	919901	361	330-495
Panel 1	MF-040518-01	167	144-215
Panel 2	MF-040518-02	52.6	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)	5847
---	------

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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 08 Jul 2020
Signature/Date