

Product Number:	101672
Lot Number:	502805
Expiration:	27-Sep-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	104012
Detector Reagent	101673	106817
SBG Reagent	101675	032904
Sample Diluent	101676	036318
Calibrator Diluent	101677	031414
Calibrator Concentrate	101734	927002
RGP Reagent	103159	N/A <sup>1</sup>

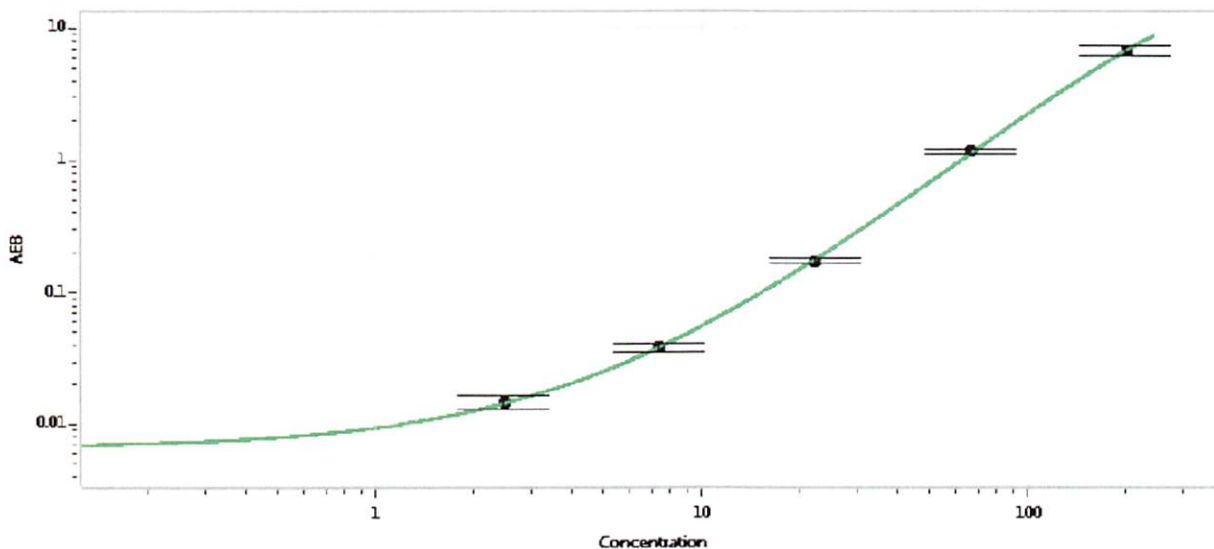
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918301	19.0	9.47-23.7
Control 2	919901	410	325-487
Panel 1	024003	127	93.6-149
Panel 2	024003	265	198-326

\* 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 18 Mar 2021  
Signature/Date