

Product Number:	101672
Lot Number:	502514
Expiration:	22-Apr-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	024006
Detector Reagent	101673	022303
SBG Reagent	101675	022305
Sample Diluent	101676	015423
Calibrator Diluent	101677	020422
Calibrator Concentrate	101734	911207
RGP Reagent	103159	N/A <sup>1</sup>

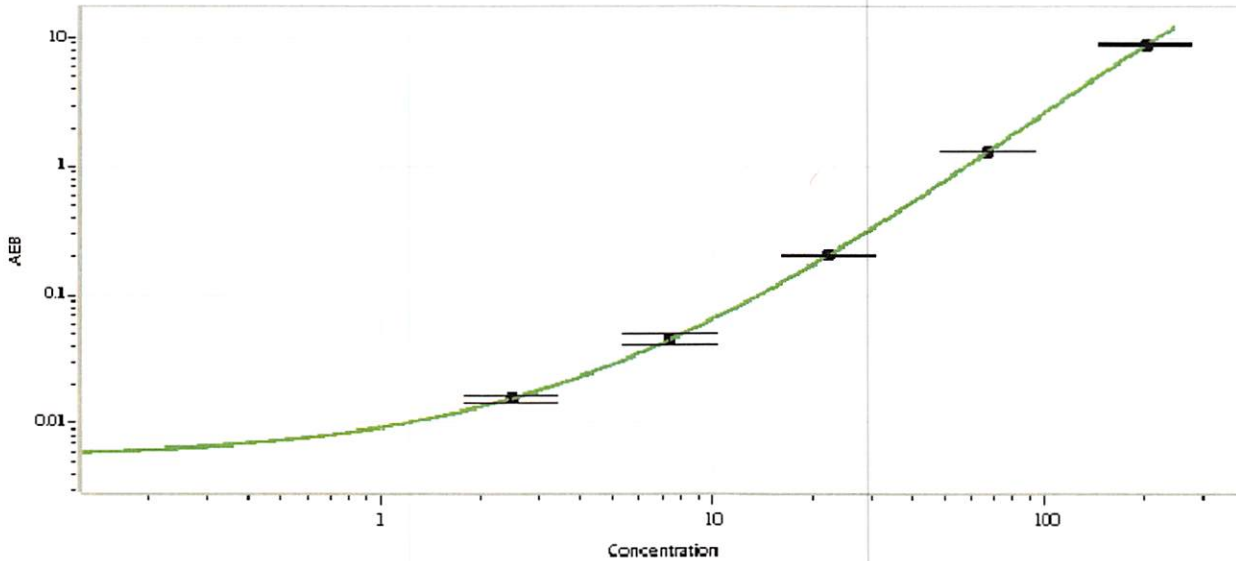
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	918301	17.1	10.8-20.7
Control 2	919901	392	330-495
Panel 1	024003	115	100-151
Panel 2	024003	255	219-329

\* 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	QC Manager	Shivani Goel 02Dec2020
	Name	Title	Signature/Date