

Product Number:	101672
Lot Number:	503150
Expiration:	24-Feb-2022
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	126803
Detector Reagent	101673	112719
SBG Reagent	101675	108710
Sample Diluent	101676	105519
Calibrator Diluent	101677	108125
Calibrator Concentrate	101734	107906
RGP Reagent	103159	N/A <sup>1</sup>

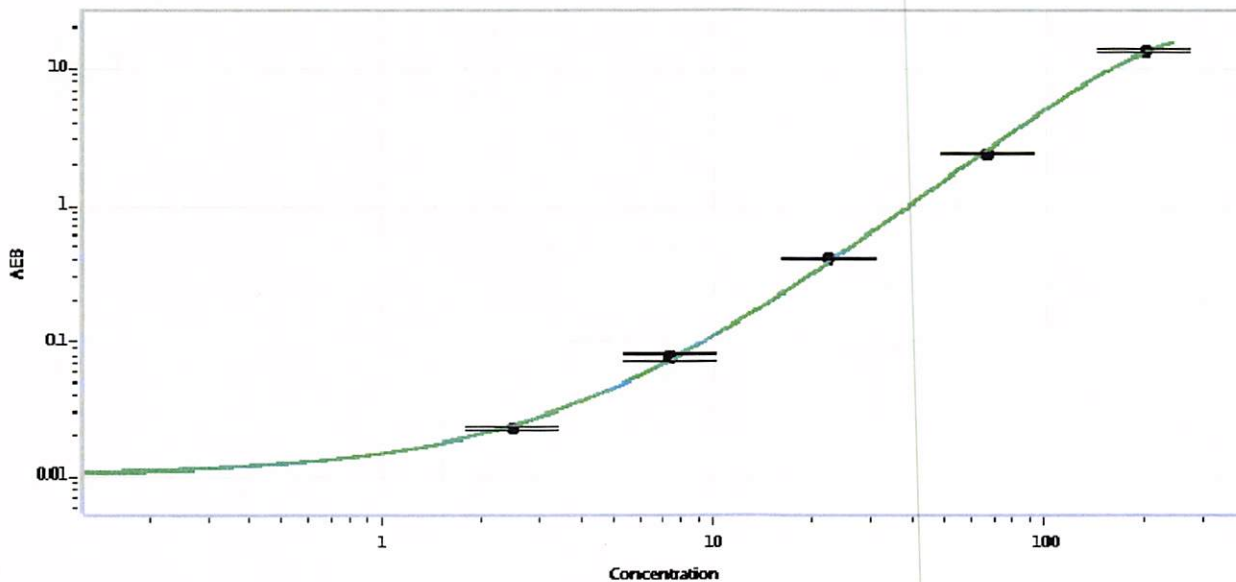
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	126904	20.4	13.2 - 23.9
Control 2	126904	368	308 - 484
Panel 1	024003	116	93.6 - 149
Panel 2	024003	244	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)	4973
---	------

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	Junpil Park	QC Supervisor		01 Nov 2021
	Name	Title	Signature/Date	