

Product Number:	101672
Lot Number:	503151
Expiration:	22-Mar-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	110804
SBG Reagent	101675	110805
Sample Diluent	101676	109024
Calibrator Diluent	101677	108131
Calibrator Concentrate	101734	107906
RGP Reagent	103159	N/A ¹

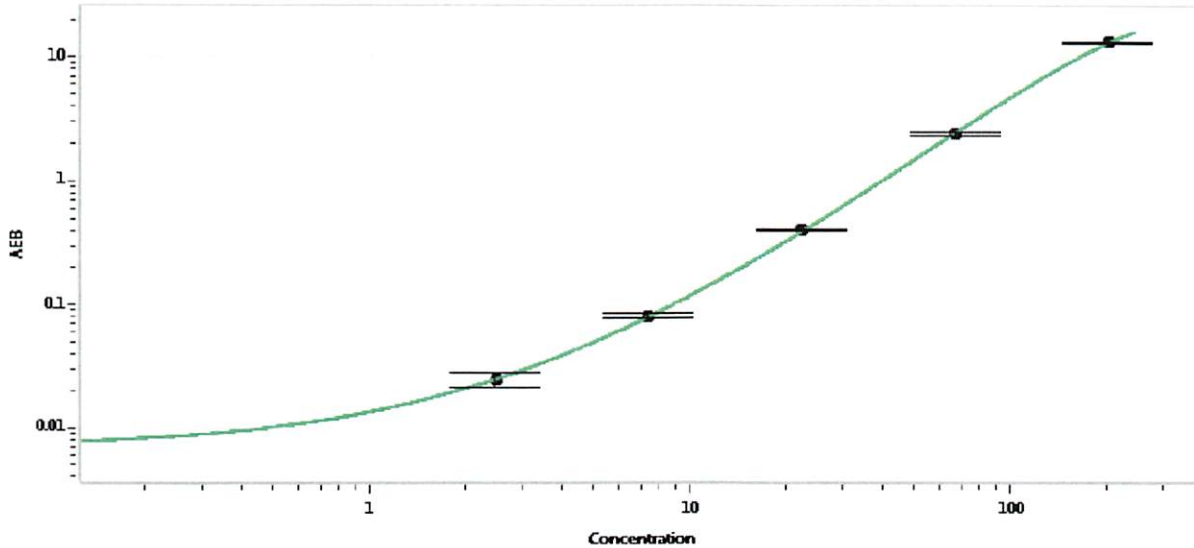
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	126904	21.6	12.5 - 23.9
Control 2	126904	435	330 - 494
Panel 1	024003	132	93.6 - 149
Panel 2	024003	270	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
Name

QC Supervisor
Title

 10 Oct 2021
Signature/Date