

Product Number:	101672
Lot Number:	503194
Expiration:	28-Jun-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	129704
Detector Reagent	101673	129803
SBG Reagent	101675	129907
Sample Diluent	101676	117927
Calibrator Diluent	101677	128016
Calibrator Concentrate	101734	130203
RGP Reagent	103159	N/A ¹

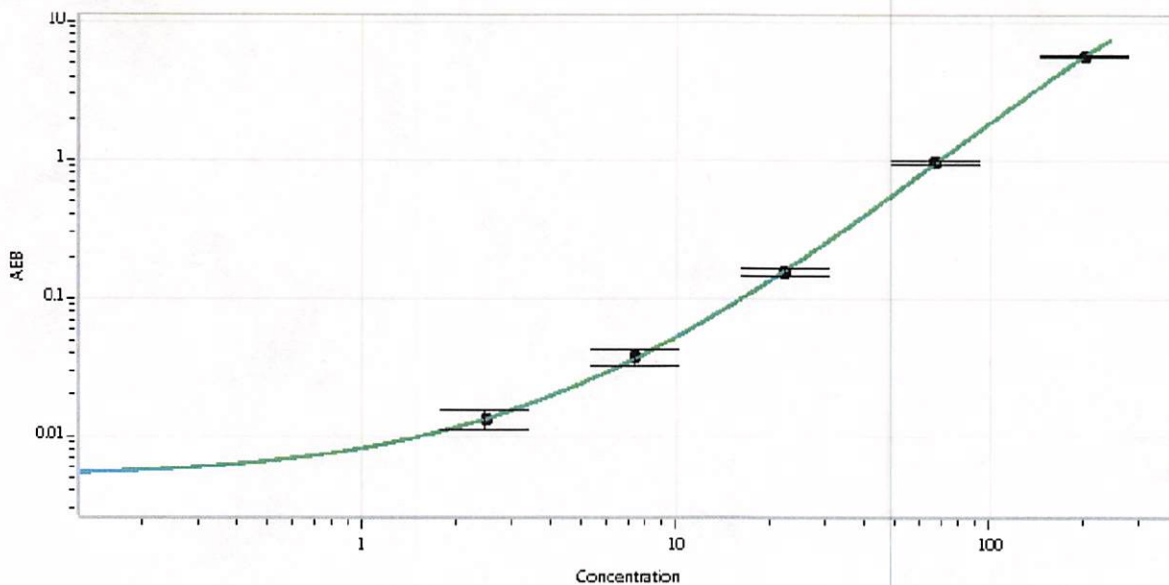
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	126904	19.2	13.2-23.9
Control 2	126904	399	308-484
Panel 1	024003	118	93.6 - 149
Panel 2	024003	261	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)	4295
---	------

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), [Signature] 03Dec2021 (Signature/Date)