

Product Number:	101132
Lot Number:	503188
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	101119	128901
Detector Reagent	101121	129002
SBG Reagent	101131	129003
Sample Diluent	103319	117926
Calibrator Diluent	101125	128014
Calibrator Concentrate	101123	129006
RGP Reagent	103159	N/A ¹

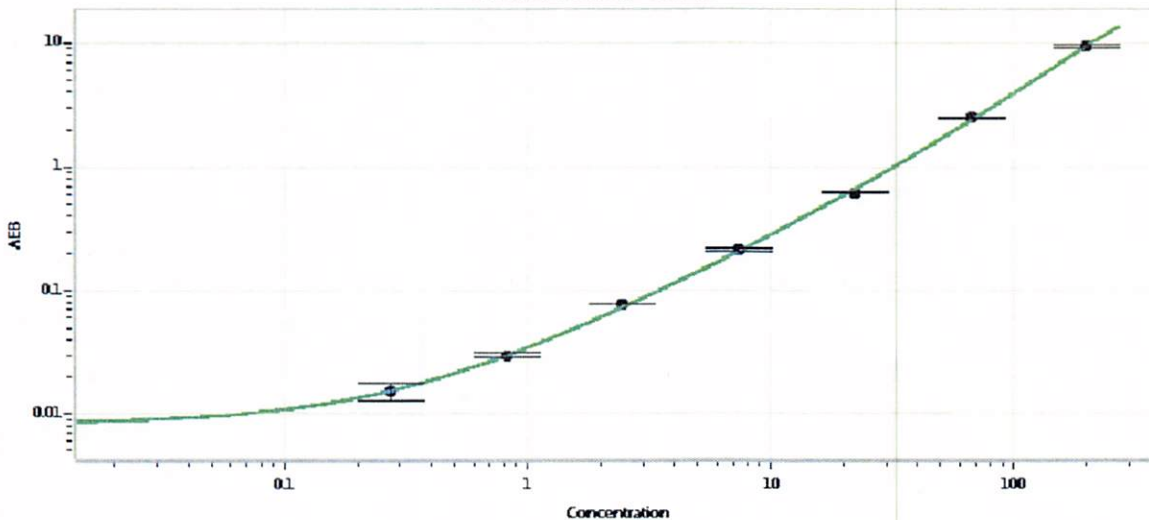
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	130713	8.41	6.70-10.1
Control 2	130713	270	154-301
Panel 1	106201	150	99.6-215
Panel 2	106201	95.4	69.9-118
Panel 3	106201	309	224-336
Panel 4	106201	259	198-296

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	6925
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval	Junpil Park	QC Supervisor		24 NOV 2021
	Name	Title	Signature/Date	