

Product Number:	101605
Lot Number:	502633
Expiration:	14-Nov-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	101607	032510
Detector Reagent	101606	032909
SBG Reagent	101608	032604
Sample Diluent	103200	032801
Calibrator Diluent	101610	031904
Calibrator Concentrate	100895	010901
RGP Reagent	103159	N/A ¹

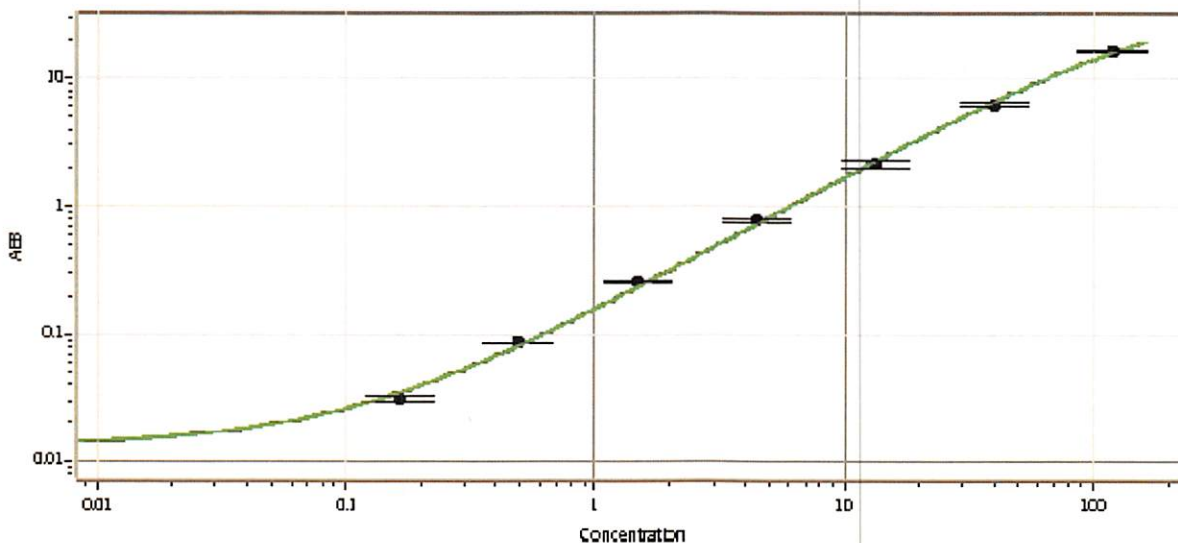
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911406	1.89	1.09 - 3.60
Control 2	911406	103	87.5 - 204
Panel 1	033901	2.83	2.13 - 3.97
Panel 2	033901	0.326	0.279 - 0.521
Panel 3	033901	19.7	19.6 - 29.4
Panel 4	033901	52.2	52.6 - 78.9

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

Calibrator Concentrate Value Assignment (pg/mL)	4866
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Junpil Park

QC Supervisor

[Signature] 15 Dec 2020

Name

Title

Signature/Date