

Product Number:	101605
Lot Number:	503096
Expiration:	2-Aug-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	101607	121902
Detector Reagent	101606	122705
SBG Reagent	101608	122104
Sample Diluent	103200	122110
Calibrator Diluent	101610	121413
Calibrator Concentrate	100895	122105
RGP Reagent	103159	N/A <sup>1</sup>

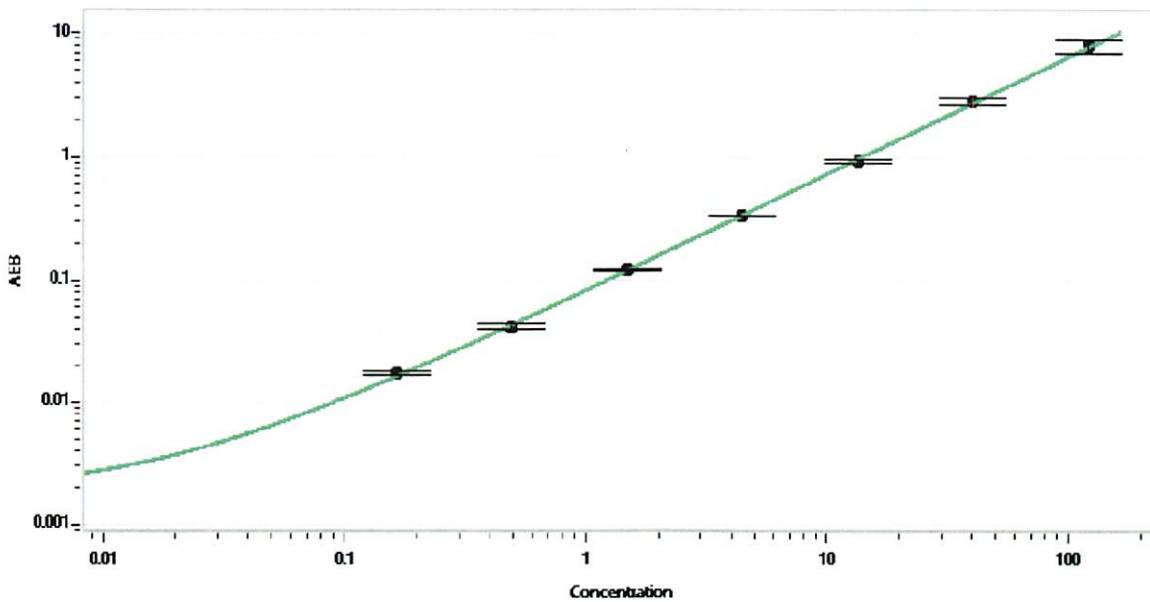
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	117610	141	90.0 - 150
Control 2	117610	2.20	1.50 - 2.50
Panel 2	033901	0.363	0.252 - 0.617
Panel 3	033901	24.0	19.1 - 31.1
Panel 4	033901	60.2	54.0 - 83.1

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

Calibrator Concentrate Value Assignment (pg/mL)	3515
---	------

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  
Name

QC Supervisor  
Title

03sep2021  
Date

Signature/Date