

Product Number:	100196
Lot Number:	502949
Expiration:	30-Oct-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	100391	112509
Detector Reagent	100315	104506
SBG Reagent	100316	104505
Sample Diluent	100317	030406
Calibrator Diluent	100665	103213
Calibrator Concentrate	100664	005203
RGP Reagent	103159	N/A ¹

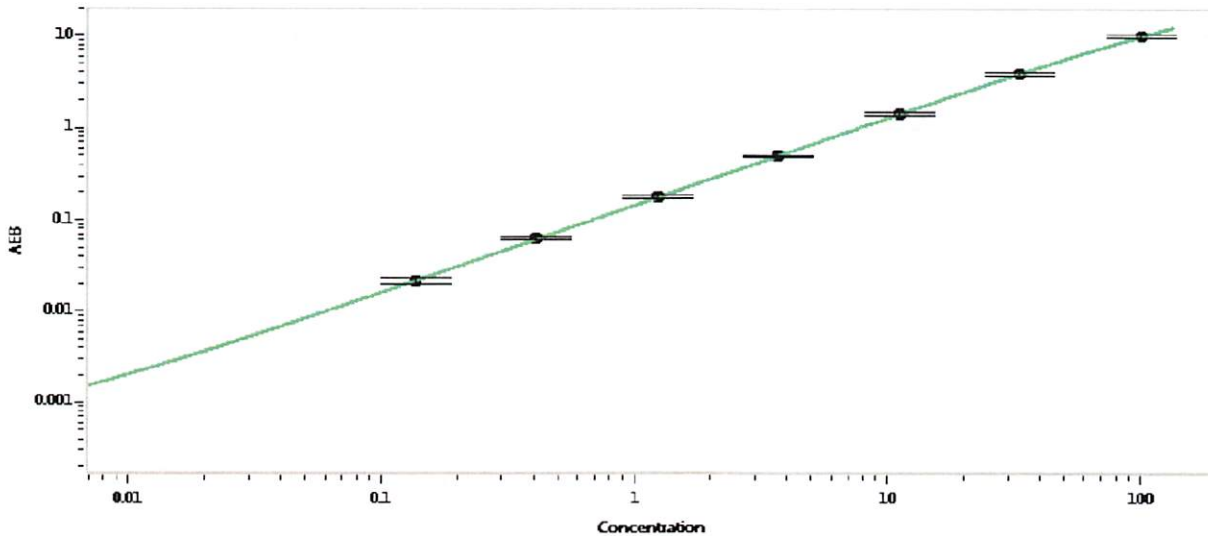
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	005203	97.5	60.0 - 100
Control 2	005203	9.50	6.00 - 10.0
Panel 1a	029002	1.54	1.17 - 2.69
Panel 2	029002	7.81	5.32 - 12.1
Panel 3	029002	81.6	49.8 - 116

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

Calibrator Concentrate Value Assignment (pg/mL)	1747
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 09 Jun 2021
Signature/Date