

Product Number:	100196
Lot Number:	503102
Expiration:	1-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	100391	122503
Detector Reagent	100315	122604
SBG Reagent	100316	122605
Sample Diluent	100317	115229
Calibrator Diluent	100665	118510
Calibrator Concentrate	100664	121803
RGP Reagent	103159	N/A ¹

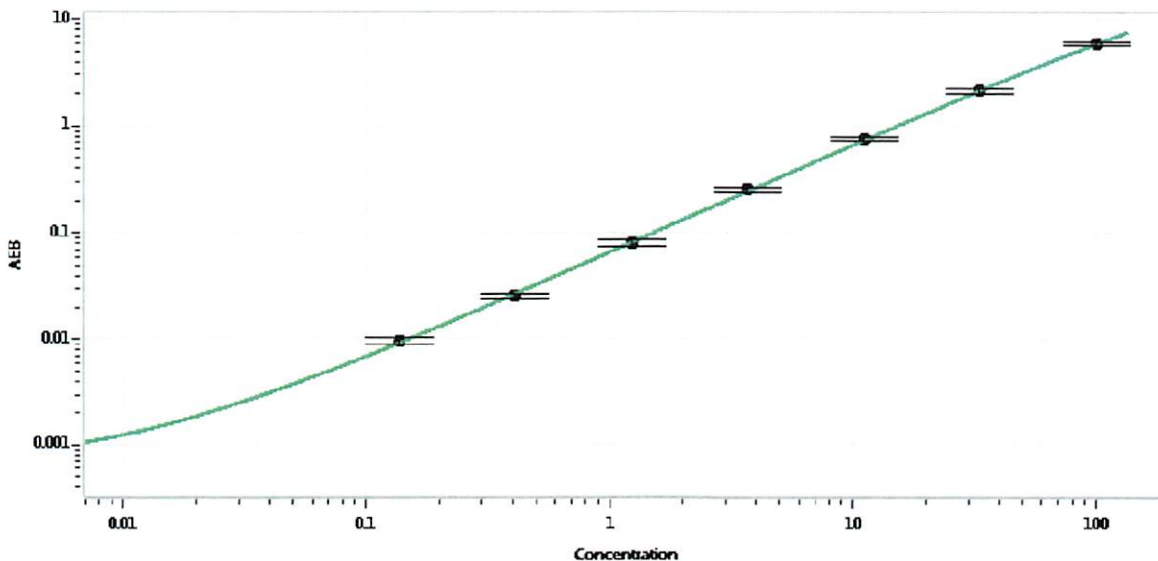
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	123410	9.61	6.41 - 10.5
Control 2	123410	97.3	65.4 - 108
Panel 1a	029002	2.28	1.17 - 2.69
Panel 2	029002	11.0	5.32 - 12.1
Panel 3	029002	85.0	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)	2009
---	------

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/Date

07Sep21