

Product Number:	100196
Lot Number:	502561
Expiration:	2-Jun-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	021903
Detector Reagent	100315	021902
SBG Reagent	100316	027604
Sample Diluent	100317	015430
Calibrator Diluent	100665	027612
Calibrator Concentrate	100664	005203
RGP Reagent	103159	N/A ¹

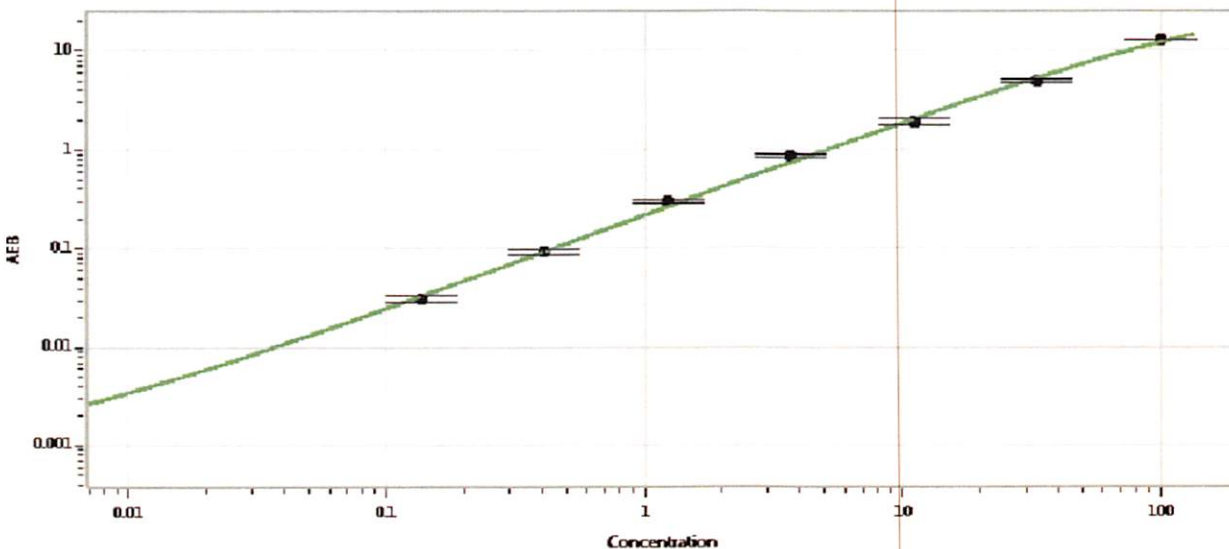
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910002	8.70	7.24-10.9
Control 2	910002	90.4	73.9-111
Panel 1a	029002	1.97	1.33-2.68
Panel 2	029002	9.34	5.33-12.4
Panel 3	029002	89.8	47.4-118

* 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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 02 NOV 2020
Signature/Date