

Product Number:	101552
Lot Number:	503137
Expiration:	18-Feb-2022
Platform(s):	HD-1/HD-x & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101554	108819
Detector Reagent	101553	125603
SBG Reagent	101555	108815
Sample Diluent	101556	104919
Calibrators	101559	112004
Control 1	101560	112005
Control 2	101561	112005
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

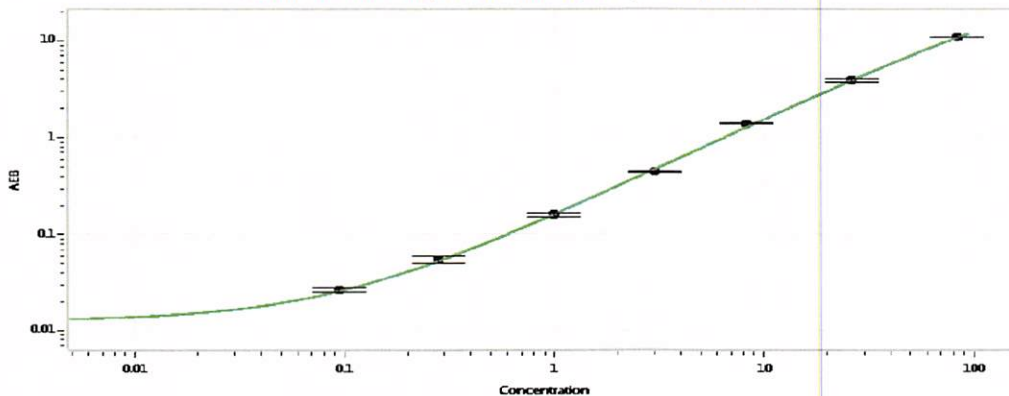
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112005	1.29	0.576 - 1.48
Control 2	112005	59.1	45.5 - 74.0
Panel 2	117205	2.98	1.90 - 5.09
Panel 3	117205	21.9	17.7 - 26.5
Panel 4	119612	35.3	21.9 - 37.9

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

Calibration Curve

2021-09-14 Tau 2.0 (Kit Lot 503137) QC (CU-81) Curve 1



Calibrator Levels (pg/mL)

A	0.000
B	0.095
C	0.281
D	1.00
E	3.01
F	8.30
G	26.0
H	83.2

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date 17 Sep 2021