

Product Number:	101599
Lot Number:	503015
Expiration:	14-Jan-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	101601	117608
Detector Reagent	102012	117804
SBG Reagent	101602	117809
Sample Diluent	103886	117810
Calibrator Diluent	102014	117116
Calibrator Concentrate	100778	001414
RGP Reagent	103159	N/A ¹

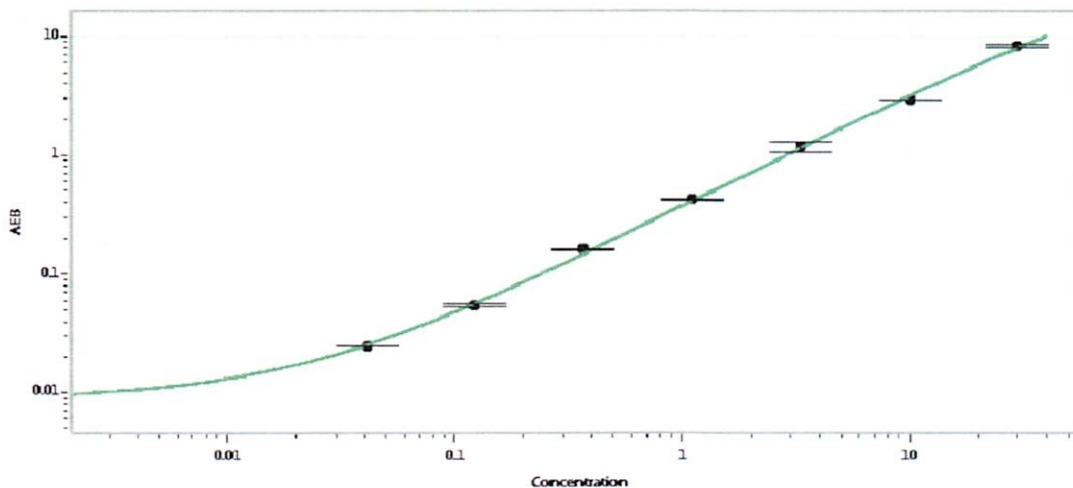
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108214	0.234	0.146-0.396
Control 2	109503	44.3	28.7-47.3
Panel 3	034908	10.9	8.12-12.5
Panel 4	034908	40.2	22.4-47.5

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

Calibrator Concentrate Value Assignment (pg/mL)	706
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Junpil Park

QC Supervisor

02 Jul 2021

Name

Title

Signature/Date