

Product Number:	101599
Lot Number:	504025
Expiration:	13-Oct-2024
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101601	328606
Detector Reagent	102012	330318
SBG Reagent	101602	329102
Sample Diluent	103886	329213
Calibrator Diluent	102014	328616
Calibrator Concentrate	100778	329107
RGP Reagent	103159	N/A ¹

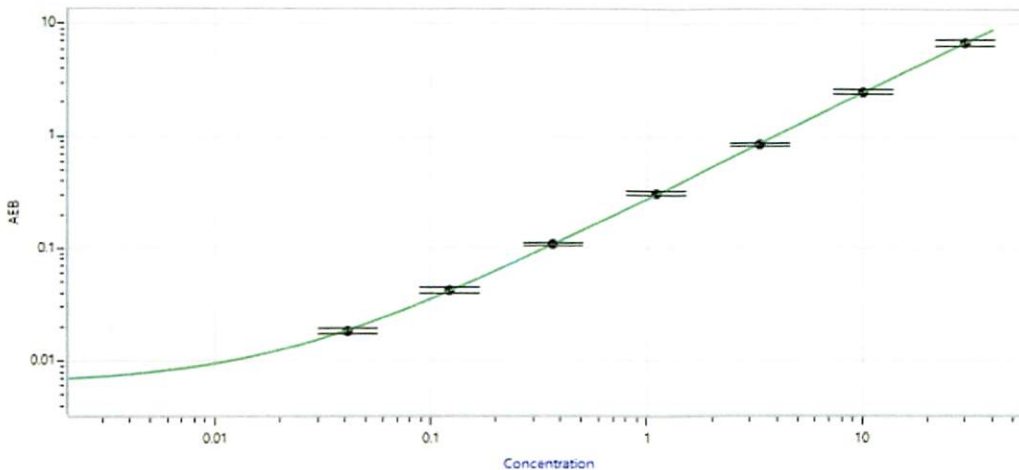
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	209305	0.446	0.432 - 0.749
Control 2	209305	26.6	9.37 - 50.1
Panel 1	321205	1.80	1.50 - 2.25
Panel 2	321205	0.308	0.250 - 0.440
Panel 3	319303	8.25	5.13 - 10.7
Panel 4	319303	29.1	20.9 - 34.1

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

Calibrator Concentrate Value Assignment (pg/mL)	418
---	-----

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

Sonia Khillan

Name

QC

Title

Sonia Khillan 06DEC2023

Signature/Date