

Product Number:	103464
Lot Number:	503889
Expiration:	24-May-2024
Platform(s):	HD-1 / HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103452	314404	314404A
Detector Reagent	103454	314503	314503A
SBG Reagent	103456	314504	314504A
Sample Diluent Reagent	103458	314605	314605A
Calibrator Diluent	103460	314606	314606A
Calibrator Concentrate	103459	316307	316307A
RGP Reagent	103159	N/A ¹	N/A ¹

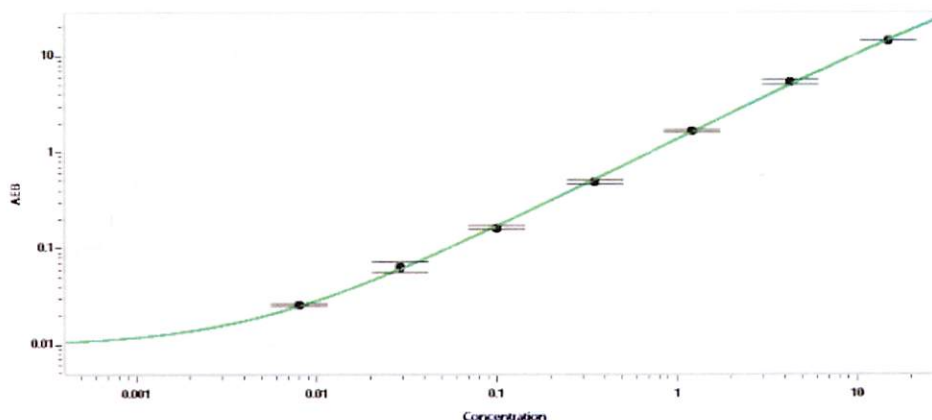
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	027502	0.166	0.134 - 0.249
Control 2	027502	10.3	7.77 - 11.7
Panel 1	106413	0.065	0.050 - 0.148
Panel 2	106413	0.030	0.019 - 0.076
Panel 3	106413	4.42	3.40 - 5.79
Panel 4	106413	17.5	10.6 - 24.9

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

Calibrator Concentrate Value Assignment (pg/mL)	345
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.008
C	0.029
D	0.100
E	0.350
F	1.22
G	4.29
H	15.0

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia Khillan 12 JUL 2023
Signature/Date