

Product Number:	103337
Lot Number:	503673
Expiration:	29-Jul-2023
Platform(s):	HD-1/HD-X & SR-X



Component	Item Number	Lot Number
Bead Reagent	103334	306116
Detector Reagent	101091	222405
SBG Reagent	101101	222406
Sample Diluent	103336	222407
Calibrator Diluent	101095	221015
Calibrator Concentrate	101093	208601
RGP Reagent	103159	N/A ¹

Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

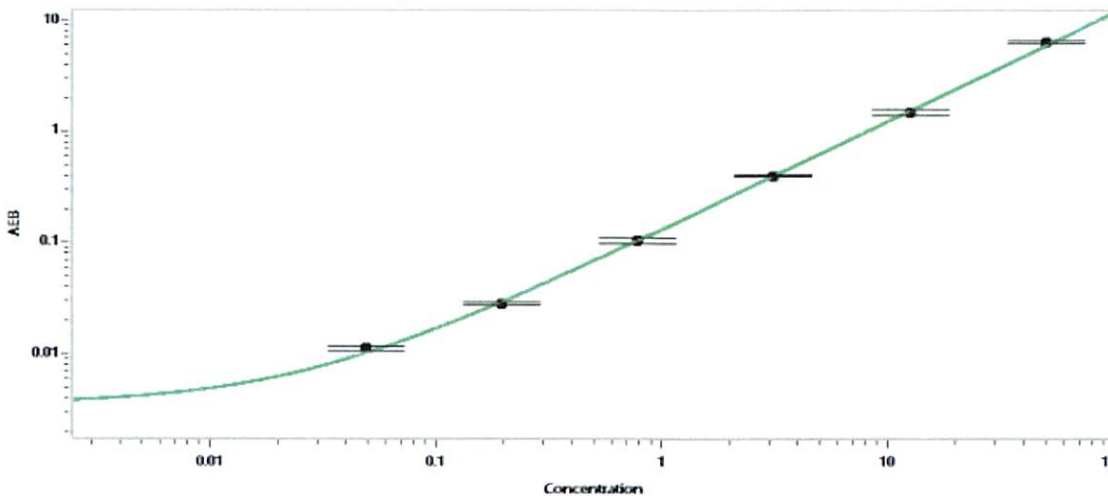
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	206704	1.54	1.13 - 1.72
Control 2	206704	31.5	21.3 - 34.0
Panel 1	211906	2.17	1.26 - 2.41
Panel 2	211906	4.11	2.37 - 4.51
Panel 3	211906	41.1	21.0 - 45.2

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

Calibrator Concentrate Value Assignment (pg/mL)	2,026
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.195
D	0.781
E	3.13
F	12.5
G	50.0

Review/Approval	<u>Sonia Khillan</u>	QC	<u>R. Della</u>	15MAR2023
	Name	Title	Signature/Date	