

Product Number:	103337
Lot Number:	503842
Expiration:	1-Feb-2024
Platform(s):	HD-1/HD-X & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	103334	306116
Detector Reagent	101091	307404
SBG Reagent	101101	307405
Sample Diluent	103336	307415
Calibrator Diluent	101095	303219
Calibrator Concentrate	101093	313003 ¹
RGP Reagent	103159	N/A ²

¹ Calibrator Concentrate Control Lot Number - 313003A

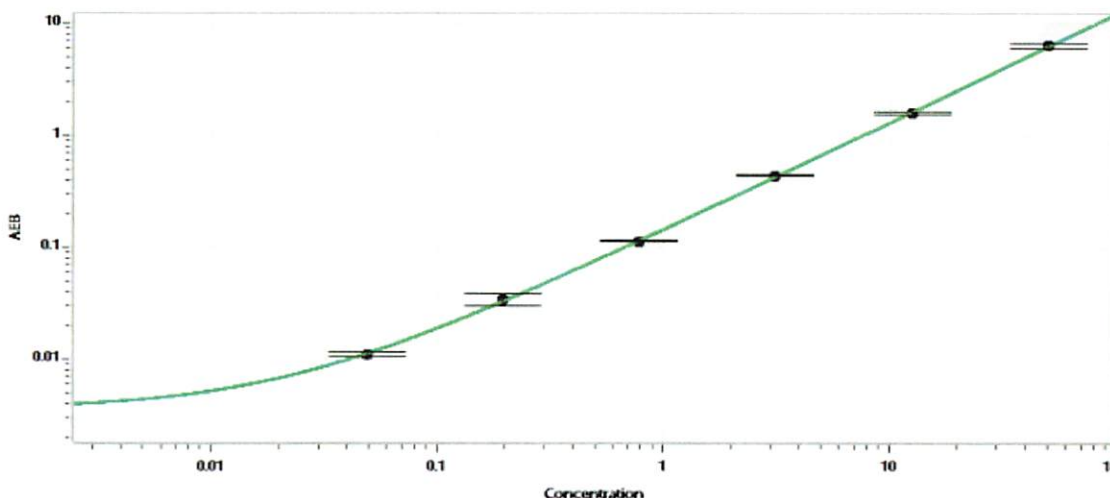
² RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	206704	1.55	1.13 - 1.72
Control 2	206704	28.4	21.3 - 34.0
Panel 1	211906	2.01	1.26 - 2.41
Panel 2	211906	3.78	2.37 - 4.51
Panel 3	211906	36.7	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)	4,352
---	-------

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: Sonia Khillan Name, QC Title, Sonia Khillan 06 JUL 2023 Signature/Date