

Product Number:	103337
Lot Number:	503741
Expiration:	18-Aug-2023
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	224208
Detector Reagent	101091	223008
SBG Reagent	101101	224507
Sample Diluent	103336	232017
Calibrator Diluent	101095	223011
Calibrator Concentrate	101093	224508
RGP Reagent	103159	N/A ¹

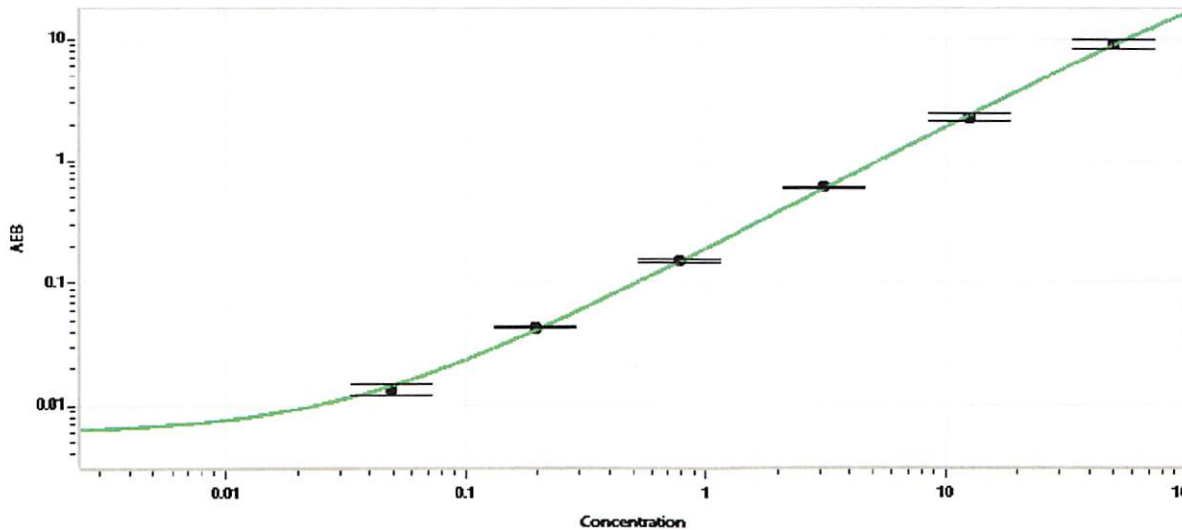
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	206704	1.59	1.07-1.82
Control 2	206704	26.8	21.6-33.1
Panel 1	211906	1.80	1.32-2.32
Panel 2	211906	3.34	2.41-4.40
Panel 3	211906	32.8	21.9-43.1

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

Calibrator Concentrate Value Assignment (pg/mL)	1,400 pg/mL
---	-------------

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 Manager Title, Sonia Khillan / 21NOV2022 Signature/Date