

Product Number:	101599
Lot Number:	503653
Expiration:	11-Apr-2023
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	213305
Detector Reagent	102012	213309
SBG Reagent	101602	213310
Sample Diluent	103886	213704
Calibrator Diluent	102014	210116
Calibrator Concentrate	100778	220304
RGP Reagent	103159	N/A ¹

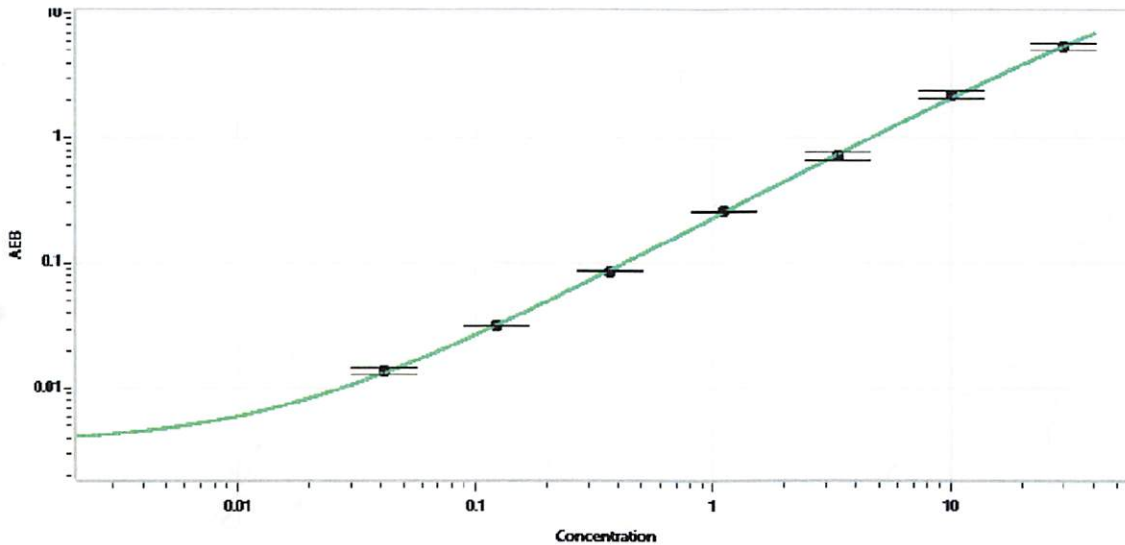
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	209305	0.549	0.411-0.729
Control 2	209305	31.1	8.21-44.0
Panel 3	034908	10.3	8.12-12.5
Panel 4	034908	34.9	22.4-47.5

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

Calibrator Concentrate Value Assignment (pg/mL)	620
---	-----

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, AC Manager Title, Sonia Khillan / 24062022 Signature/Date