

Product Number:	101599
Lot Number:	503502
Expiration:	22-May-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	216203
Detector Reagent	102012	220101
SBG Reagent	101602	216302
Sample Diluent	103886	216504
Calibrator Diluent	102014	214209
Calibrator Concentrate	100778	213307
RGP Reagent	103159	N/A ¹

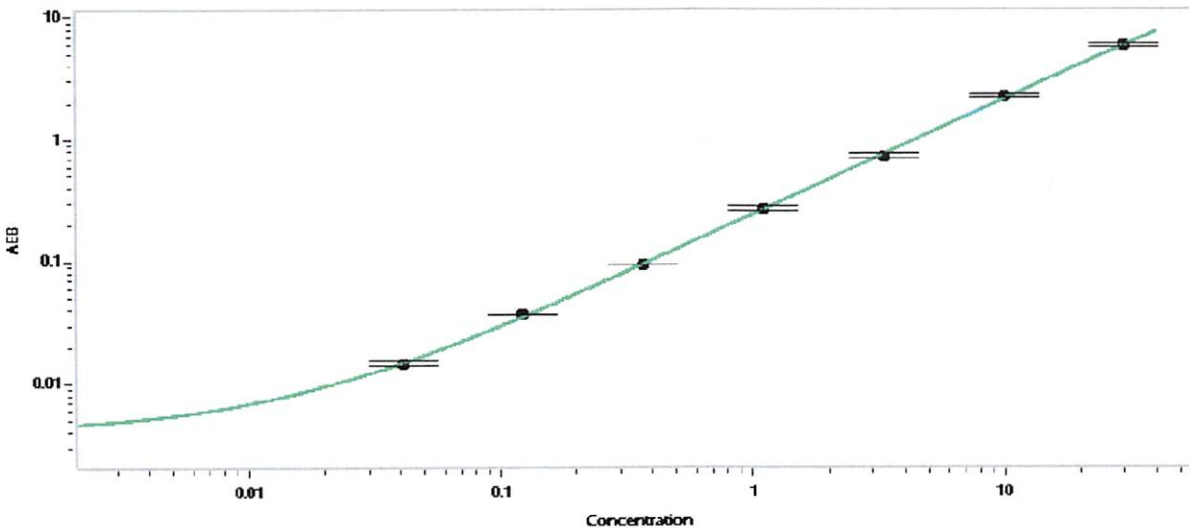
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	213307	23.4	18.0-30.0
Control 2	213307	0.449	0.375-0.625
Panel 3	034908	10.1	8.12-12.5
Panel 4	034908	36.7	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)	757
---	-----

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 QC Manager Sonia Khillan / 17AUG2022
 Name Title Signature/Date