

Product Number:	101599
Lot Number:	503736
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	231805
SBG Reagent	101602	216302
Sample Diluent	103886	216504
Calibrator Diluent	102014	214209
Calibrator Concentrate	100778	213307
RGP Reagent	103159	N/A <sup>1</sup>

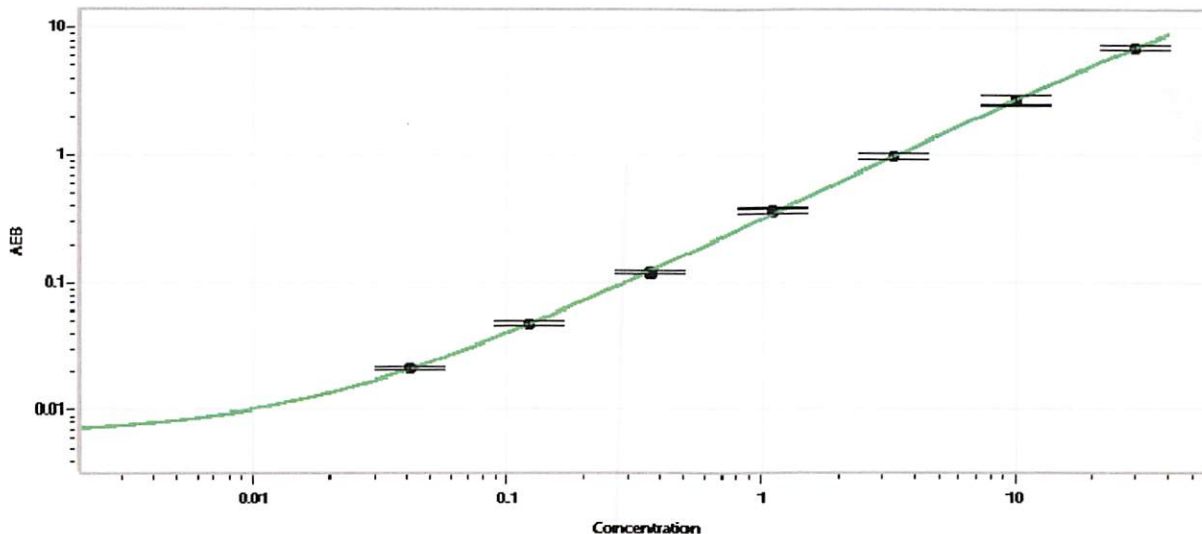
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	209305	0.647	0.418-0.752
Control 2	209305	40.8	7.13-51.0
Panel 3	034908	11.3	8.12-12.5
Panel 4	034908	43.0	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 pg/mL
---	-----------

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

QC  
Title

*Sonia Khillan* 30 NOV 2022  
Signature/Date