



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

Product Number:	101599
Lot Number:	503903
Expiration:	15-May-2024
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number	Control Number
Bead Reagent	101601	316410	316410A
Detector Reagent	102012	316309	316309A
SBG Reagent	101602	317314	317314A
Sample Diluent	103886	316404	316404A
Calibrator Diluent	102014	313510	313510A
Calibrator Concentrate	100778	319813	319813A
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>

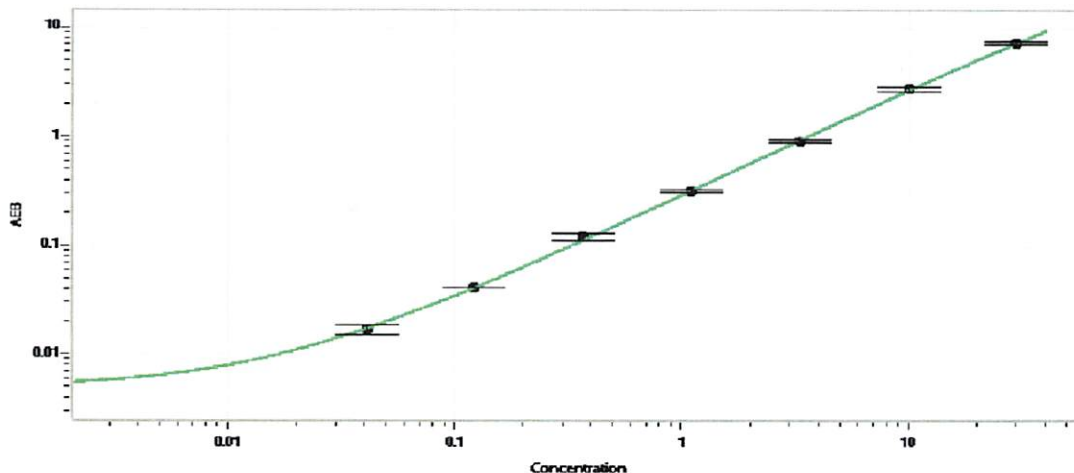
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	209305	0.616	0.418 - 0.752
Control 2	209305	33.2	7.13 - 51.0
Panel 1	321205	1.80	1.52 - 2.29
Panel 2	321205	0.333	0.271 - 0.444
Panel 3	319303	9.52	4.82 - 10.8
Panel 4	319303	31.2	20.1 - 34.0

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

Calibrator Concentrate Value Assignment (pg/mL)	506
---	-----

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

*ROBERT SOMOGYI*  
Name

QC, MANAGER  
Title

*[Signature]* 07 AUG 2023  
Signature/Date