

Product Number:	100988
Lot Number:	503829
Expiration:	13-Feb-2024
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	100974	318004	318004A
Detector Reagent	100976	305508	N/A
SBG Reagent	100987	305509	N/A
Sample Diluent	103734	305903	N/A
Calibrator Diluent	100980	304414	N/A
Calibrator Concentrate	100978	319309	319309A
RGP Reagent	103159	N/A ¹	N/A ¹

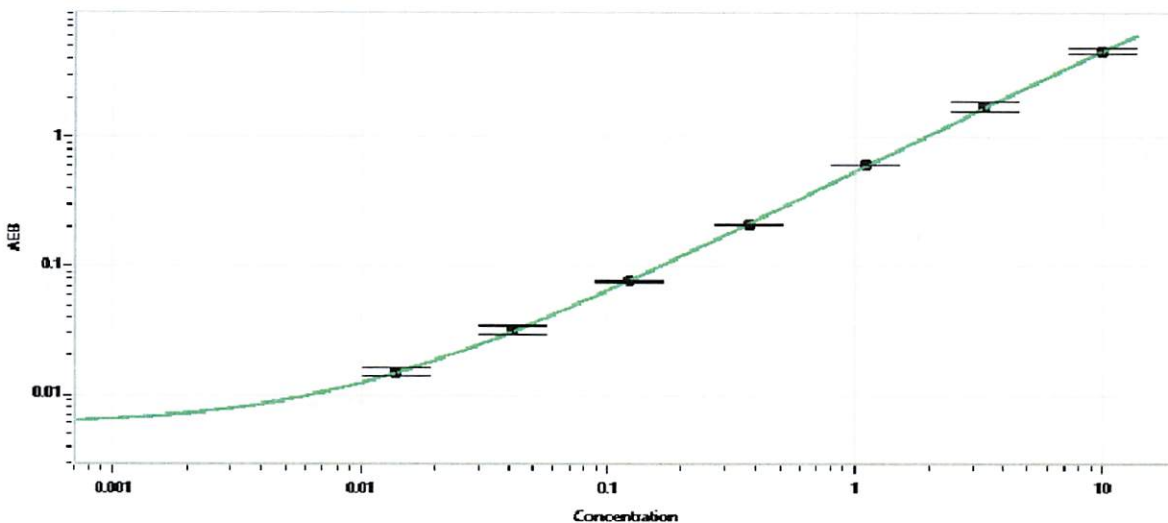
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	108314	0.149	0.079 - 0.247
Control 2	108314	13.0	10.3 - 15.7
Panel 1	030001	0.163	0.105 - 0.234
Panel 3	030001	29.5	19.9 - 38.7
Panel 4	030001	0.165	0.111 - 0.215

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

Calibrator Concentrate Value Assignment (pg/mL)	194
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

[Signature] 25JUL2023

Signature/Date