

Product Number:	101605
Lot Number:	503806
Expiration:	10-Aug-2023
Platform(s):	HD-1 / HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101607	222202
Detector Reagent	101606	222408
SBG Reagent	101608	222409
Sample Diluent	103200	222410
Calibrator Diluent	101610	303712
Calibrator Concentrate	100895	222412
RGP Reagent	103159	N/A ¹

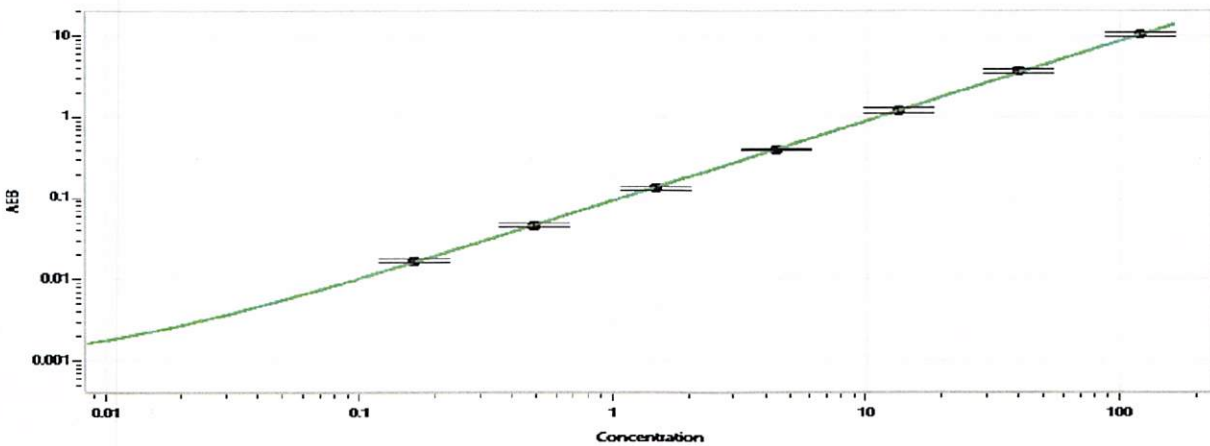
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	222412	2.31	1.50 - 2.50
Control 2	222412	141.0	90.0 - 150
Panel 1	124703	1.06	0.870 - 1.68
Panel 2	124703	1.19	0.902 - 1.94
Panel 3	124703	8.79	6.42 - 13.0
Panel 4	124703	47.8	35.8 - 77.4

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

Calibrator Concentrate Value Assignment (pg/mL)	2596
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Matt Koulobonis

QC supervisor

M. Koulobonis 09Feb2023

Name

Title

Signature/Date

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

techsupport@quantix.com

©Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures