

Product Number:	101580
Lot Number:	503933
Expiration:	14-Jul-2024
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104236	319504	319504A
Detector Reagent	104327	319805	319805A
SBG Reagent	104329	319806	319806A
Sample Diluent	104331	319506	319506A
Calibrator Diluent	101585	319804	319804A
Calibrator Concentrate	104332	320906	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

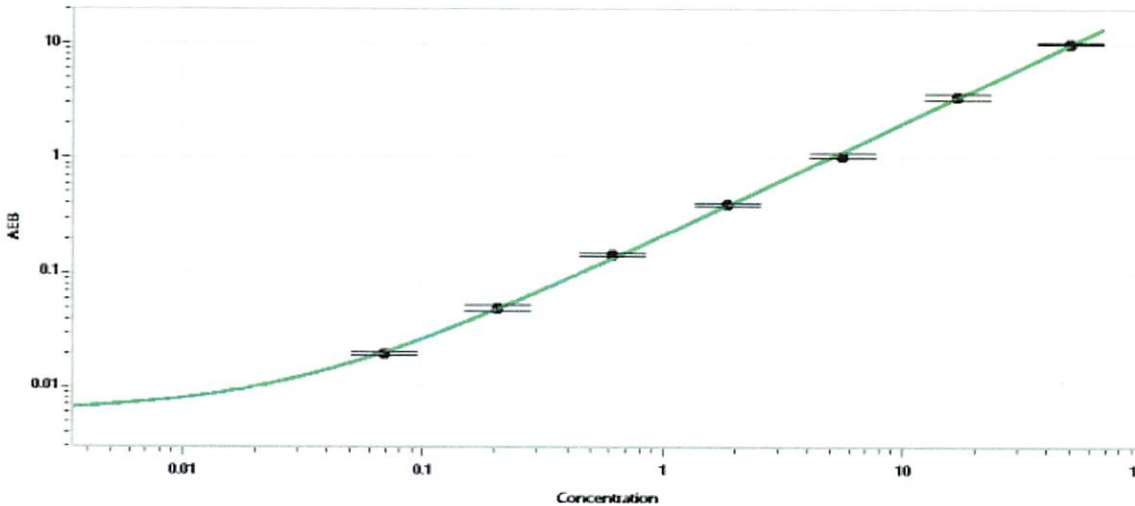
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	206502	3.20	2.21 - 3.51
Control 2	206502	52.4	35.6 - 58.8
Panel 2	115317	2.32	1.51 - 3.73
Panel 3	116301	30.6	15.5 - 39.3

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

Calibrator Concentrate Value Assignment (pg/mL)	1408
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval


QC
 17 AUG 2023

Name
Title
Signature/Date