

Product Number:	104111
Lot Number:	503907
Expiration:	1-Aug-2024
Platform(s):	HD-1 / HD-X & SR-X



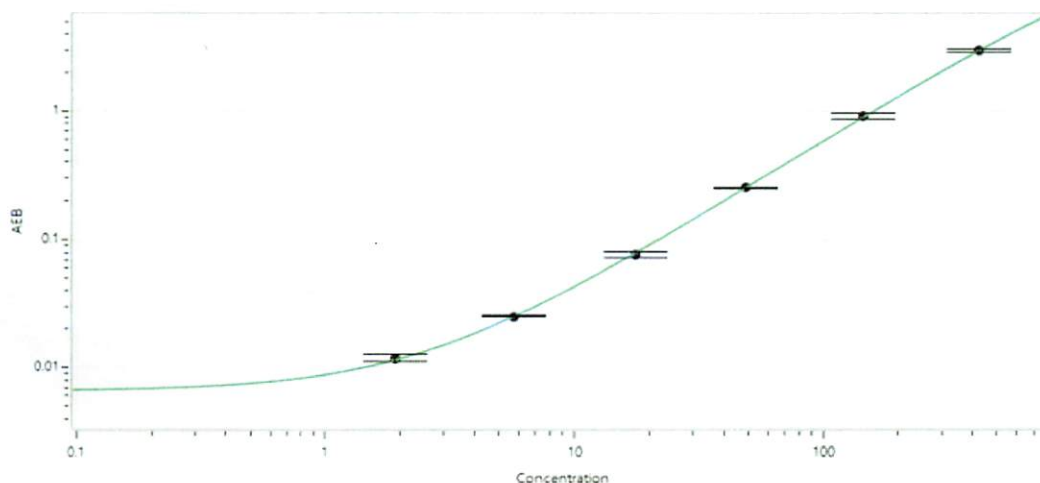
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	104098	321404
Detector Reagent	104099	322205
SBG Reagent	104100	321306
Sample Diluent	104101	321309
Reference Calibrator	104106	322604
Control 1	104107	322605
Control 2	104108	322605
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1	322605	42.6	34.0 - 51.0	
Control 2	322605	725	614 - 922	
Panel 1	301702	16.5	12.7 - 20.9	
Panel 2	301702	14.6	12.3 - 20.9	
Panel 3	309704	452	370 - 661	
Panel 4	301702	96.5	57.9 - 125	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.91
C	5.70
D	17.7
E	49.1
F	144
G	419

Review/Approval

Sonia Khillgn
Name

QC

Title

Sumit K. 20SEP2023
Signature/Date