

Product Number:	104111
Lot Number:	503843
Expiration:	10-Apr-2024
Platform(s):	HD-1 / HD-X & SR-X



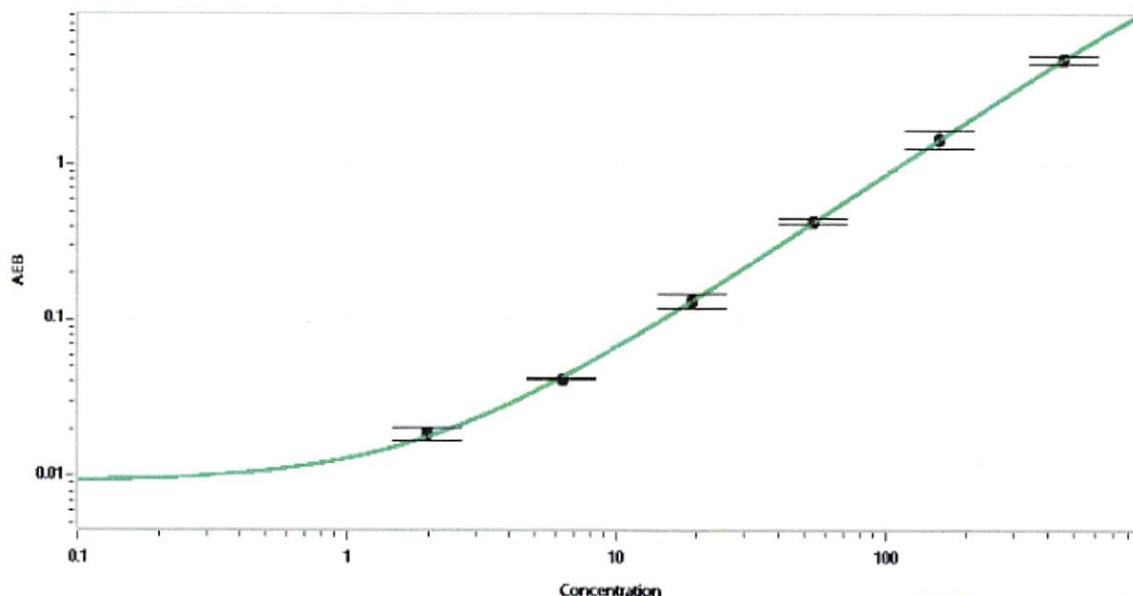
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104098	310012	N/A
Detector Reagent	104099	310307	310307A
SBG Reagent	104100	310103	310103A
Sample Diluent	104101	310104	310104A
Reference Calibrator	104106	311402	311402A
Control 1	104107	311403	311403A
Control 2	104108	311403	311403A
RGP Reagent	103159	N/A ¹	N/A

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	311403	46.4	38.2 - 57.3	
Control 2	311403	916	729 - 1093	
Panel 1	301702	16.1	13.4 - 20.2	
Panel 2	301702	15.2	13.2 - 20.0	
Panel 3	309704	449	429 - 644	
Panel 4	301702	94.2	64.7 - 119	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.98
C	6.28
D	19.1
E	53.6
F	158
G	454

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia Khillan 22 MAY 2023
Signature/Date