

Product Number: 100860
Lot Number: 503926
Expiration: 25-Jul-2024
Platform(s): HD-1 & HD-X



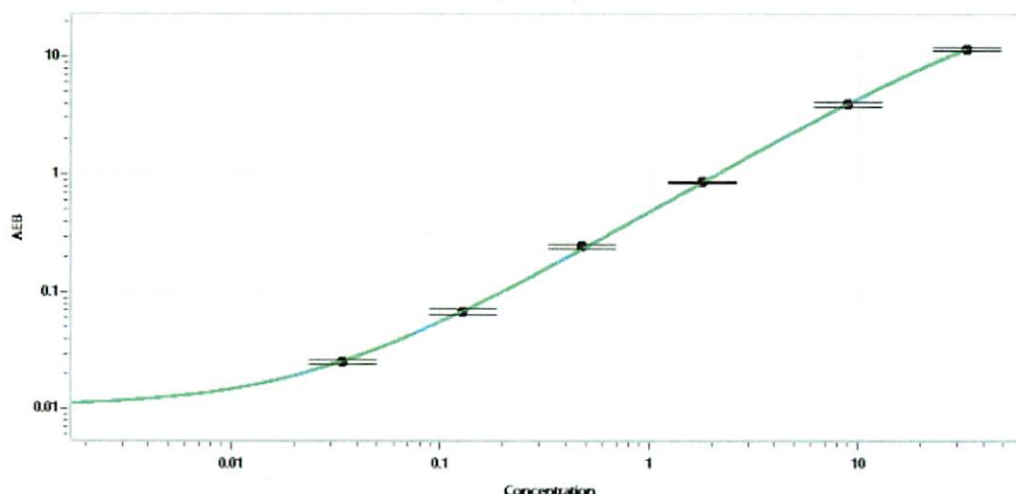
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	320710
Detector Reagent	100862	321208
SBG Reagent	100868	323008
Sample Diluent	100870	320605
Calibrators	101111	320902
Control 1	100875	320903
Control 2	100876	320903
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	320903	6.04	5.23 - 8.10	
Control 2	320903	0.097	0.075 - 0.127	
Panel 1	306704	0.794	0.772 - 1.16	
Panel 2	306704	3.94	2.98 - 7.27	
Panel 3	306704	15.7	10.1 - 25.4	

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.034
C	0.128
D	0.481
E	1.79
F	8.98
G	32.9

Review/Approval: Sonia Khillan QC Sonia Khillan 25SEP2023
Name Title Signature/Date