

Product Number:	100860
Lot Number:	503527
Expiration:	18-Jul-2023
Platform(s):	HD-1 & HD-X



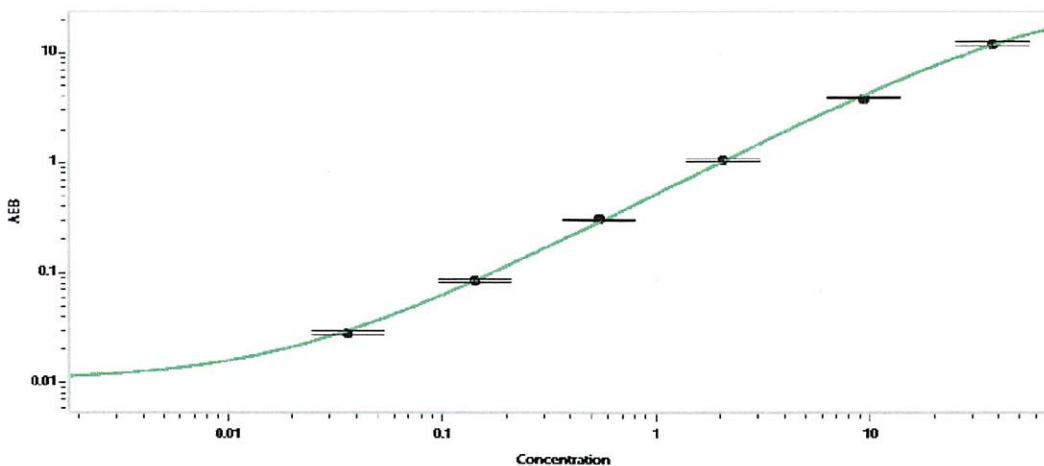
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	219908
Detector Reagent	100862	220011
SBG Reagent	100868	229104
Sample Diluent	100870	220013
Calibrators	101111	213303
Control 1	100875	213304
Control 2	100876	213304
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	213304	6.43	5.07-7.61	
Control 2	213304	0.102	0.089-0.140	
Panel 1	115916	1.23	0.830-1.35	
Panel 2	115916	6.80	1.78-9.41	
Panel 3	115916	43.5	22.1-52.6	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.036
C	0.142
D	0.537
E	2.04
F	9.35
G	37.2

Review/Approval Sonia Khillan Name QC Manager Title Sonia Khillan / 26.07.2022 Signature/Date