

Product Number:	100860
Lot Number:	503868
Expiration:	10-Apr-2024
Platform(s):	HD-1 & HD-X



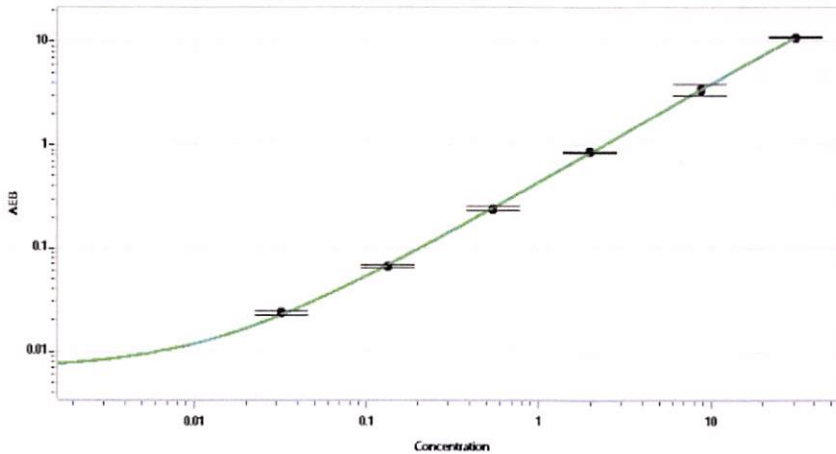
Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number	Control Number
Bead Reagent	100865	310303	310303A
Detector Reagent	100862	310710	310710A
SBG Reagent	100868	310404	310404A
Sample Diluent	100870	310005	N/A
Calibrators	101111	310701	310701A
Control 1	100875	310702	310702A
Control 2	100876	310702	310702A
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	310702	6.51	4.92 - 7.34
Control 2	310702	0.114	0.081 - 0.125
Panel 1	306704	0.863	0.772 - 1.16
Panel 2	306704	5.11	2.98 - 7.27
Panel 3	306704	17.5	8.18 - 26.3

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



Calibrator Levels (pg/mL)

A	0.000
B	0.032
C	0.132
D	0.538
E	1.98
F	8.64
G	30.5

Review/Approval Sonia Khillan QC Sonia Khillan 02MAY2023
 Name Title Signature/Date