

Product Number:	100860
Lot Number:	503376
Expiration:	4-May-2023
Platform(s):	HD-1 & HD-X



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

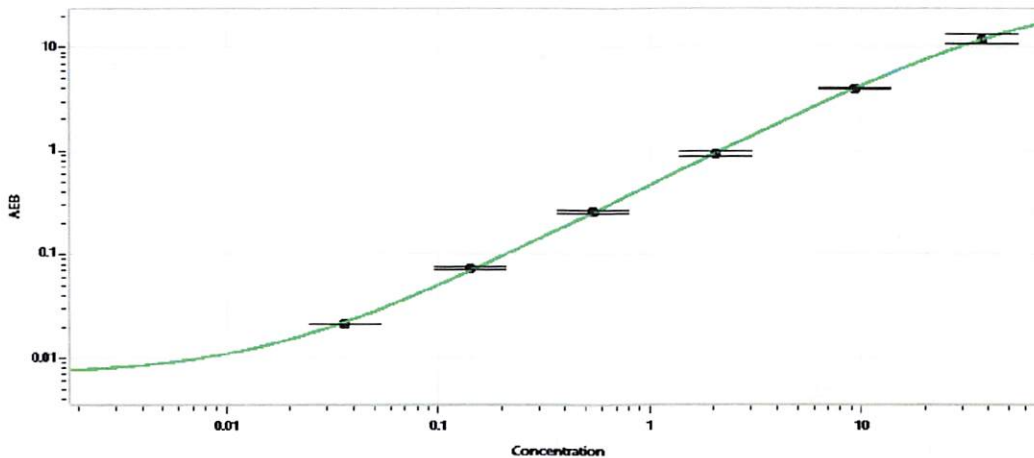
Component	Part Number	Lot Number
Bead Reagent	100865	212602
Detector Reagent	100862	212405
SBG Reagent	100868	212406
Sample Diluent	100870	212407
Calibrators	101111	213303
Control 1	100875	213304
Control 2	100876	213304
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	213304	6.50	5.05-7.58
Control 2	213304	0.127	0.094-0.141
Panel 1	115916	1.14	0.830-1.35
Panel 2	115916	4.47	1.78-9.41
Panel 3	115916	34.3	22.1-52.6

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

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 / 07SEP2022 Signature/Date