

Product Number:	100860
Lot Number:	504691
Expiration:	14-Nov-2025
Platform(s):	HD-1/HD-X



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

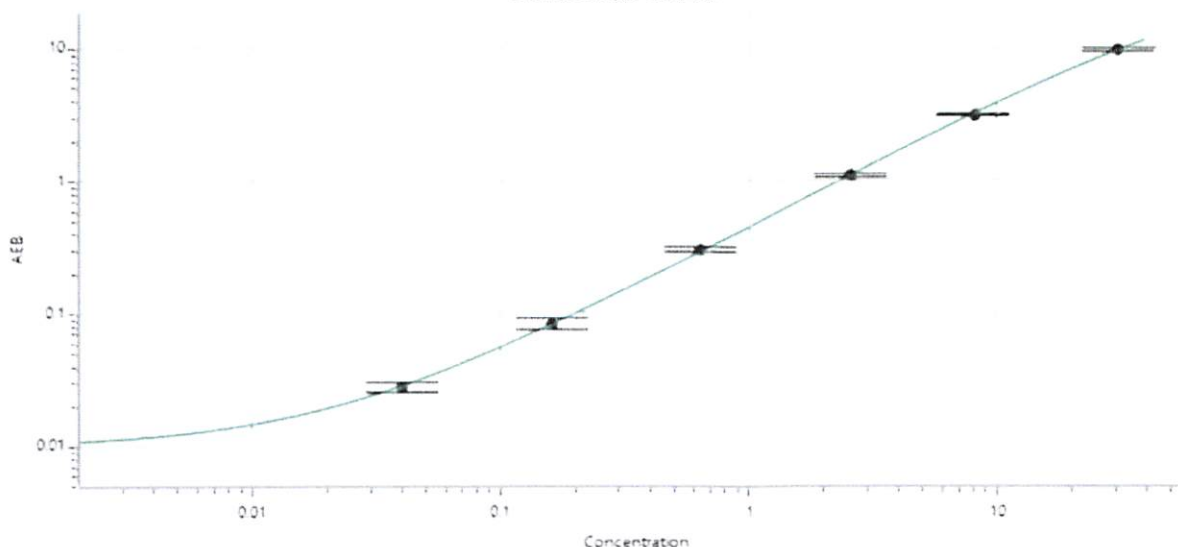
Component	Part Number	Lot Number
Bead Reagent	100865	522606
Detector Reagent	100862	522605
SBG Reagent	100868	522603
Sample Diluent	100870	522306
Reference Calibrators	101111	331808
Control 1	100875	331809
Control 2	100876	331809
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331809	5.58	4.56 - 6.84
Control 2	331809	0.120	0.096 - 0.145

* 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	0.040
C	0.161
D	0.634
E	2.55
F	8.06
G	30.5

Review/Approval

Dhruvi Patel QA Specialist II, QA

Name

Title

EDP | 27 Aug 2025

Signature/Date