

Product Number:	100860
Lot Number:	504592
Expiration:	9-Oct-2026
Platform(s):	HD-1/HD-X



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

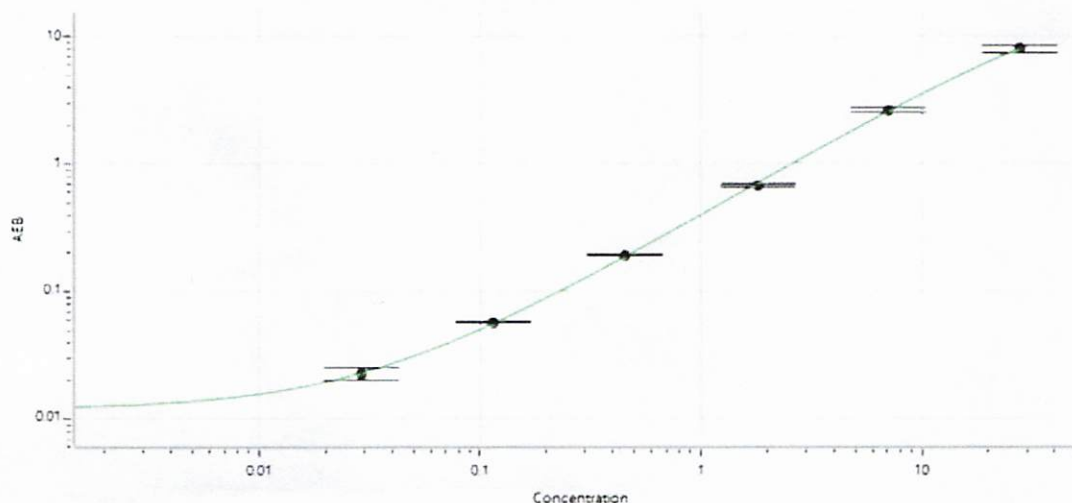
Component	Part Number	Lot Number
Bead Reagent	100865	510403
Detector Reagent	100862	510108
SBG Reagent	100868	510005
Sample Diluent	100870	509909
Reference Calibrators	101111	510506
Control 1	100875	510507
Control 2	100876	510507
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	510507	4.60	4.01 - 6.01
Control 2	510507	0.077	0.066 - 0.099

* 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	0.029
C	0.115
D	0.454
E	1.82
F	7.02
G	27.7

Review/Approval

Dilish Patel
Name

Sr. Doc. control-QA
Title

at 13 Jun 2021
Signature/Date