

Product Number:	103845	
Lot Number:	504649	
Expiration Date:	11-Dec-2025	
Platform(s):	HD-X	
Component	Part Number	Lot Number
Bead Reagent	103758	416301
Detector Reagent	103760	416511
SBG Reagent	103762	419211
Sample Diluent	103763	416302
Reference Calibrators	103766	416909
Control 1	103767	416910
Control 2	103768	416910
Control 3	103773	416910
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific



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

Release Materials*					* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (ng/mL)	Mean Low Limit (ng/mL)	Mean High Limit (ng/mL)	
Control 1	416910	306	254	459	
Control 2	416910	2671	2229	4395	
Control 3	416910	12472	9364	20443	

Review/Approval

Dhruvi Patel

Doc Control II, QA

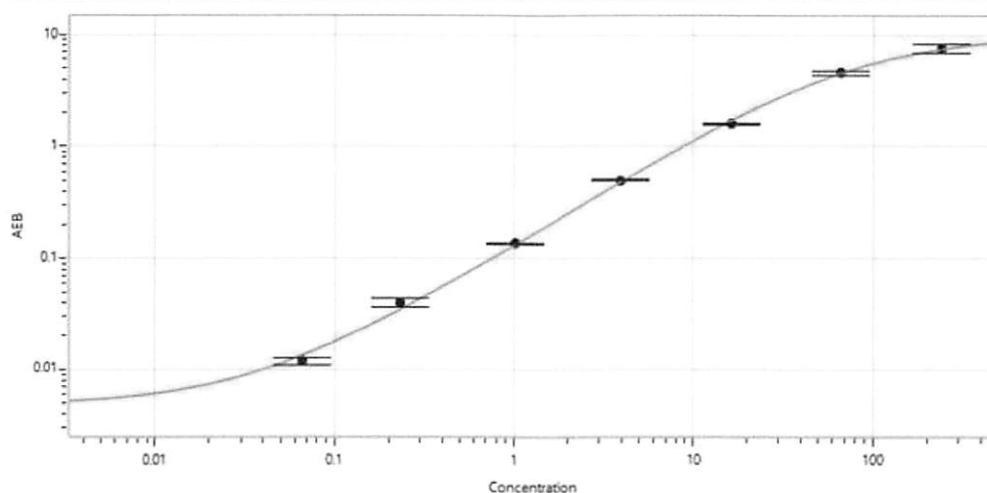
DP / 30 Jun 2025

Name

Title

Signature/Date

Calibrator Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.066
C	0.233
D	1.02
E	3.92
F	16.3
G	66.8
H	244