

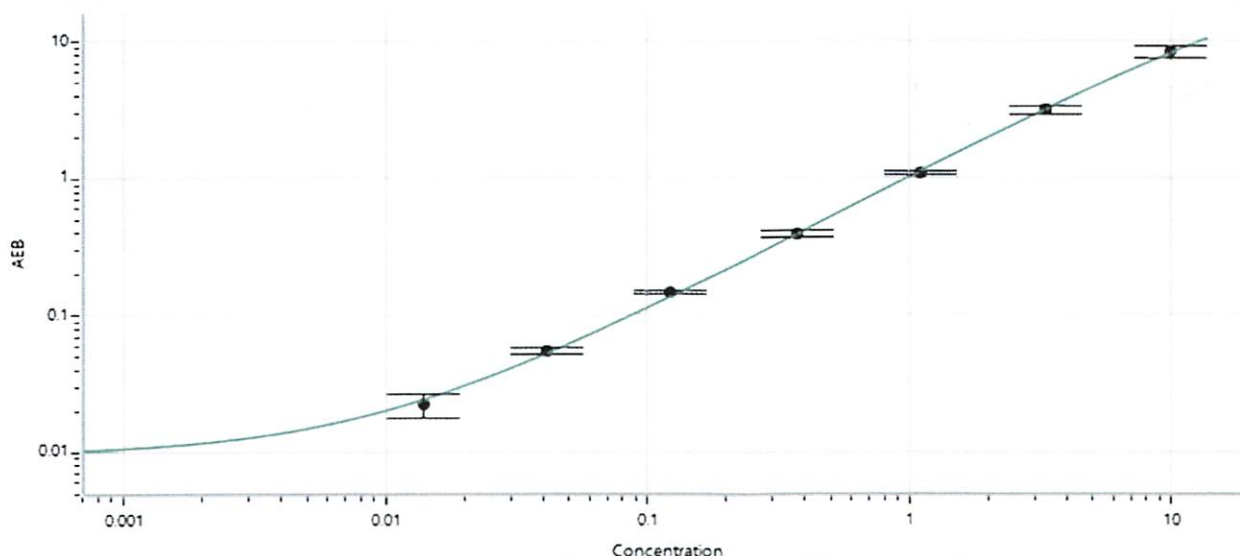
Data below represents results generated on the Simoa® HD-X Analyzer.

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403912	7.33	6.00 - 10.0
Control 2	403912	0.085	0.075 - 0.125

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

175



A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Concentration

Dipesh Patel Ser. Doc Control-QA at 09 APR 2025

Name Title Signature/Date