

Product Number:	103872
Lot Number:	504645
Expiration Date:	28-Feb-2026
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103794	516902
Detector Reagent	103796	424104
SBG Reagent	103798	424105
Sample Diluent	103800	424106
Reference Calibrators	103803	424812
Control 1	103804	424813
Control 2	103805	424813
RGP Reagent	103159	N/A ¹

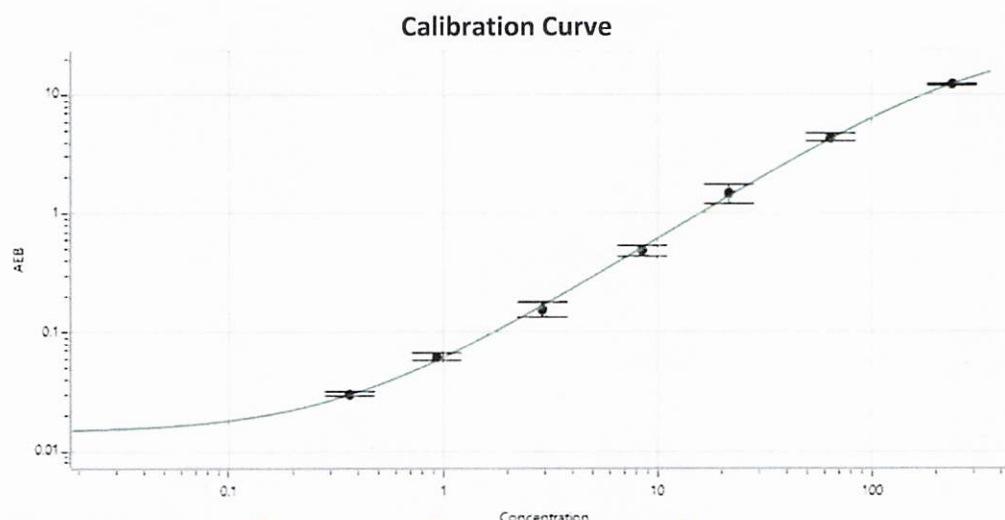
¹ Reagents are not Kit Lot Specific



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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	424813	7.35	6.51 - 9.77
Control 2	424813	148.0	125 - 187

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



Calibrator Levels (pg/mL)

A	0.000
B	0.367
C	0.927
D	2.90
E	8.49
F	21.5
G	64.4
H	236

Review/Approval

Dipesh Patel
Name

Sr. Dir. Control - QA
Title

26 Jun 2025
Signature/Date