

Product Number:	103464
Lot Number:	504609
Expiration:	5-Nov-2026
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	103452	512511
Detector Reagent	103454	513604
SBG Reagent	103456	513602
Sample Diluent	103458	512614
Calibrator Diluent	103460	512615
Calibrator Concentrate	103459	512502
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

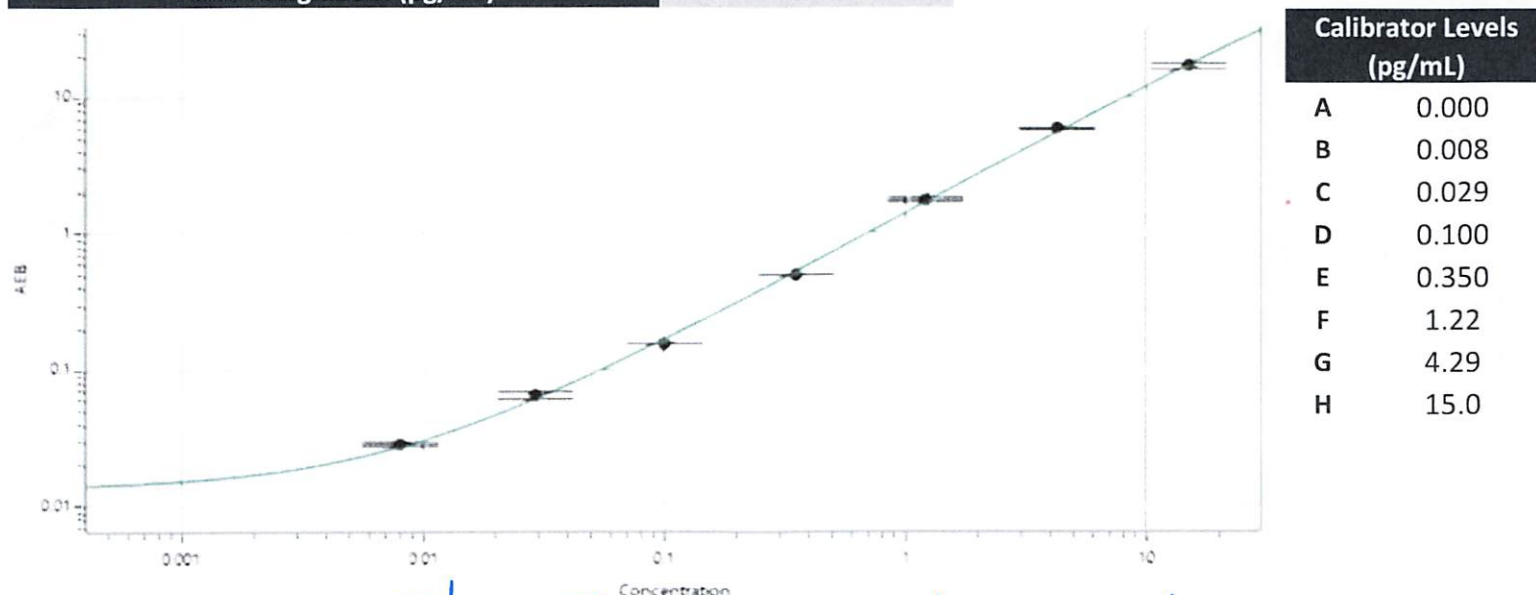


Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	027502	0.167	0.123 - 0.260
Control 2	027502	10.4	6.80 - 12.6

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

Calibrator Concentrate Value Assignment (pg/mL)	260
---	-----



Review/Approval

Dhruvi Patel

Name

DOC Control/II, QA

Title

DP / 11 Jun 2025

Signature/Date