

Product Number:	103277
Lot Number:	504690
Expiration:	5-Jan-2026
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	103265	418703
Detector Reagent	103268	418704
SBG Reagent	103269	418705
Sample Diluent	103270	418706
Calibrator Diluent	103266	418710
Calibrator Concentrate	103272	420401
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

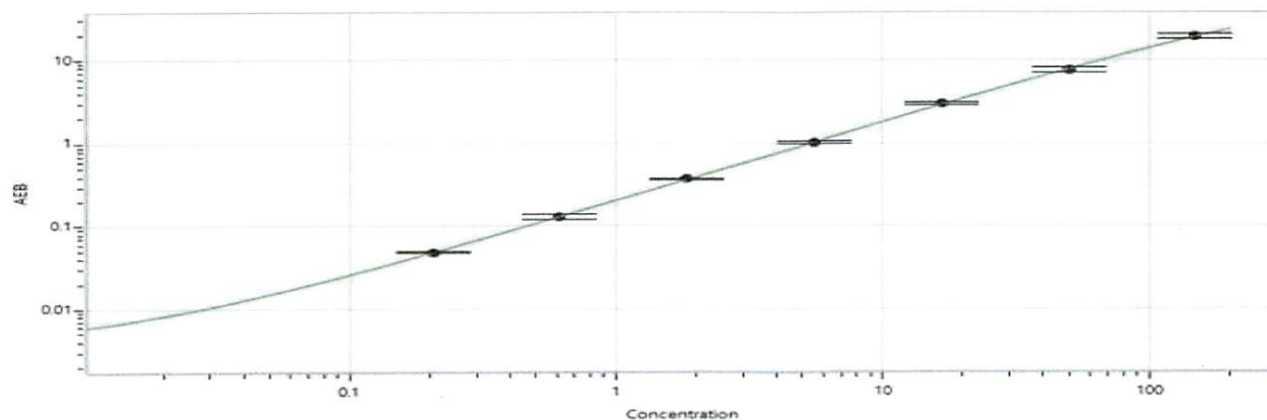


Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	420401	3.58	3.00 - 5.00
Control 2	420401	96.5	75.0 - 125

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

Calibrator Concentrate Value Assignment (pg/mL)	4933
---	------



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval

Dhruvi Patel QA Specialist II, QA *SDP* 21 AUG 2025
 Name Title Signature/Date