

Product Number:	105259
Lot Number:	504424
Expiration:	20-Feb-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	105239	429512
Detector Reagent	105241	429602
SBG Reagent	105243	429603
Sample Diluent	105244	429521
Calibrator Diluent	105246	423328
Control Diluent	105247	423327
Calibrator Concentrate	105255	411426
Control 1 Concentrate	105257	411428
Control 2 Concentrate	105258	411429
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

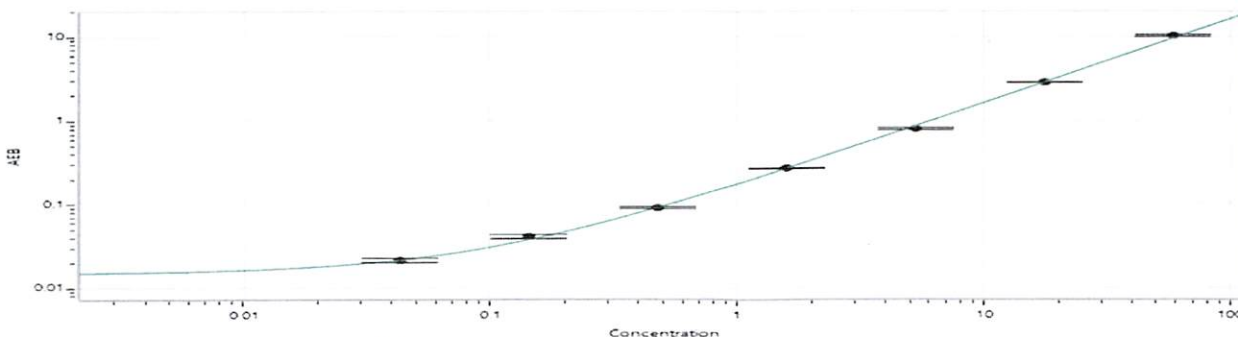
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411428	1.19	0.938 - 1.74
Control 2 Concentrate	411429	65.6	46.2 - 85.8

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

Calibrator Concentrate Reconstitution Volume (uL)	1122
Control 1 Concentrate Reconstitution Volume (uL)	1122
Control 2 Concentrate Reconstitution Volume (uL)	1122

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.043
C	0.143
D	0.477
E	1.59
F	5.29
G	17.6
H	58.7

Review/Approval

Dhruvi Patel *DOC Control I, QA* *DP* 10 Dec 2024

Name

Title

Signature/Date