

Product Number:	104997
Lot Number:	504533
Expiration:	9-Oct-2026
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104982	509904
Detector Reagent	104984	510406
SBG Reagent	104986	510407
Sample Diluent	104987	509910
Calibrator Diluent	104988	509912
Control Diluent	105402	509911
Calibrator Concentrate	104993	412338
Control 1 Concentrate	104995	412340
Control 2 Concentrate	104996	412341
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



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

Release Materials*			
IL-4®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.768	0.547 - 1.02
Control 2 Concentrate	412341	19.8	14.1 - 26.1

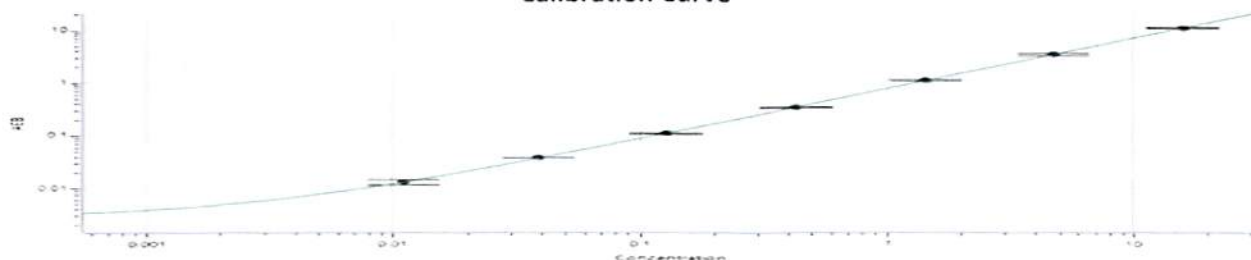
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	875
Control 1 Concentrate Reconstitution Volume (uL)	875
Control 2 Concentrate Reconstitution Volume (uL)	875

Calibrator Levels (pg/mL)

A	0.000
B	0.011
C	0.039
D	0.128
E	0.428
F	1.43
G	4.74
H	15.8

Calibration Curve



Review/Approval

Dhruvi Patel QA Specialist II, QA *DR* 19 Sep 2025
 Name Title Signature/Date

Release Materials*

IL-13®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.749	0.518 - 0.961
Control 2 Concentrate	412341	21.1	14.2 - 26.3

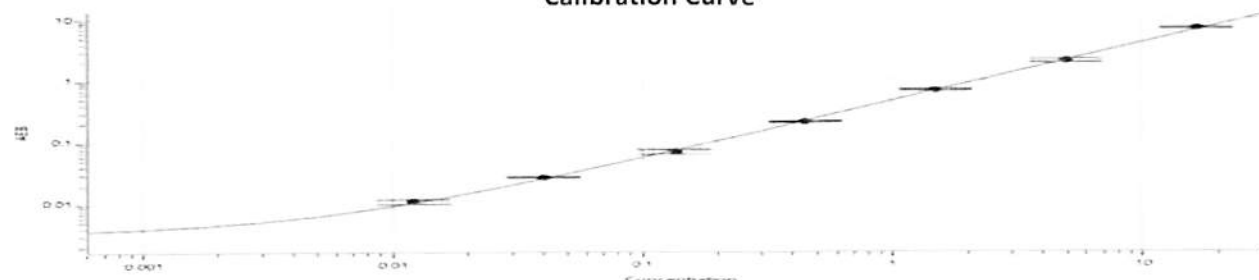
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	875
Control 1 Concentrate Reconstitution Volume (uL)	875
Control 2 Concentrate Reconstitution Volume (uL)	875

Calibrator Levels (pg/mL)

A	0.000
B	0.012
C	0.040
D	0.135
E	0.446
F	1.49
G	4.96
H	16.5

Calibration Curve



Release Materials*

IL-5®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.750	0.524 - 0.974
Control 2 Concentrate	412341	19.1	13.5 - 25.0

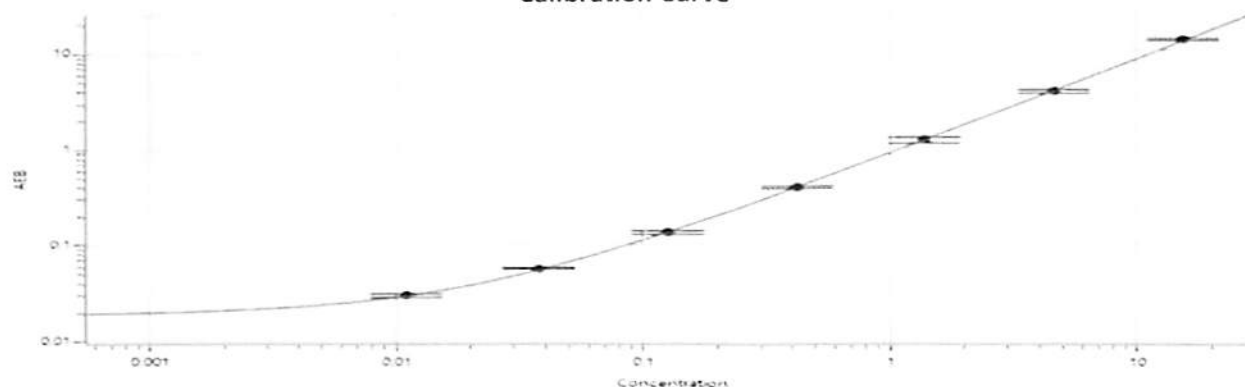
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	875
Control 1 Concentrate Reconstitution Volume (uL)	875
Control 2 Concentrate Reconstitution Volume (uL)	875

Calibrator Levels (pg/mL)

A	0.000
B	0.011
C	0.038
D	0.125
E	0.417
F	1.39
G	4.61
H	15.4

Calibration Curve



Release Materials*

IL-17A®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.569	0.395 - 0.734
Control 2 Concentrate	412341	14.3	10.0 - 18.6

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

Calibrator Concentrate Reconstitution Volume (uL)	875
Control 1 Concentrate Reconstitution Volume (uL)	875
Control 2 Concentrate Reconstitution Volume (uL)	875

**Calibrator Levels
(pg/mL)**

A	0.000
B	0.009
C	0.028
D	0.093
E	0.311
F	1.04
G	3.46
H	11.5

Calibration Curve

