

Product Number:	104997
Lot Number:	504395
Expiration:	7-Feb-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104982	425504
Detector Reagent	104984	426001
SBG Reagent	104986	426002
Sample Diluent	104987	422022
Calibrator Diluent	104988	423326
Control Diluent	105402	423325
Calibrator Concentrate	104993	412315
Control 1 Concentrate	104995	412316
Control 2 Concentrate	104996	412317
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

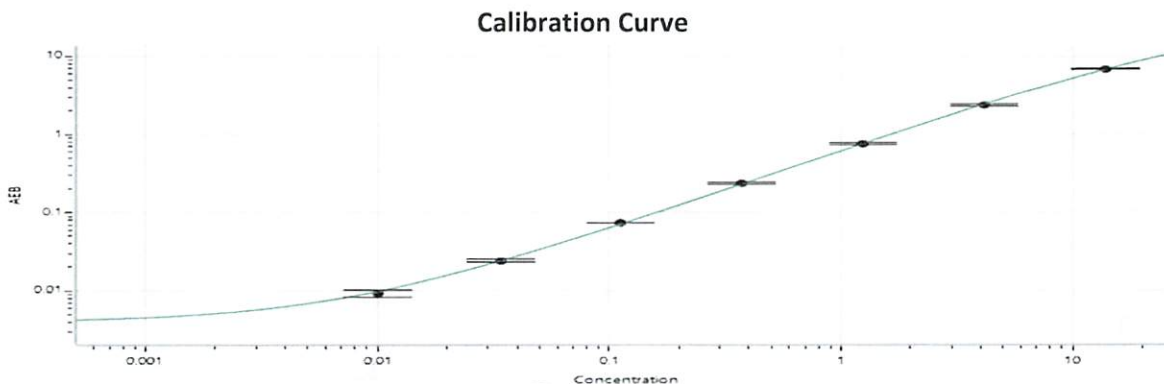
¹ Reagents are not Kit Lot Specific

Release Materials*			
IL-4®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412316	0.721	0.501 - 0.930
Control 2 Concentrate	412317	15.4	12.5 - 23.3

* 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	0.010
C	0.034
D	0.112
E	0.374
F	1.25
G	4.15
H	13.8



Review/Approval

Dipesh Patel
Name

Director of ILQA
Title

[Signature] 07 Nov 2024
Signature/Date

Release Materials*

IL-13®

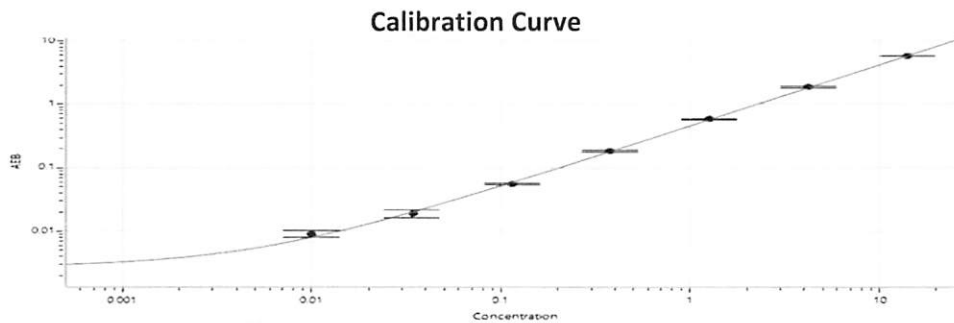
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412316	0.676	0.482 - 0.895
Control 2 Concentrate	412317	17.1	11.8 - 21.9

* 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	0.010
C	0.034
D	0.114
E	0.378
F	1.26
G	4.20
H	14.0



Release Materials*

IL-5®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412316	0.754	0.486 - 0.902
Control 2 Concentrate	412317	17.9	12.2 - 22.7

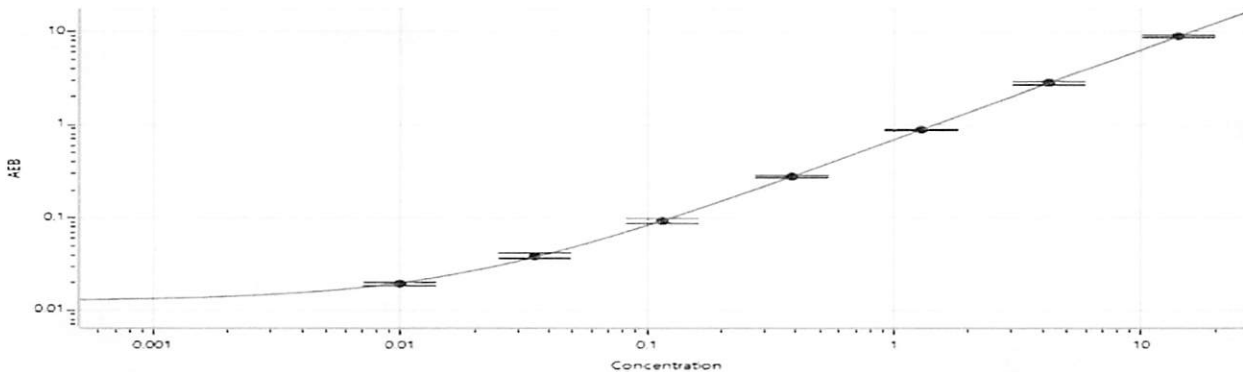
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	0.010
C	0.035
D	0.115
E	0.385
F	1.28
G	4.26
H	14.2

Calibration Curve



Release Materials*

IL-17A®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412316	0.537	0.352 - 0.653
Control 2 Concentrate	412317	11.3	8.64 - 16.0

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

Concentrate Reconstitution Volume (uL)	875
Concentrate Reconstitution Volume (uL)	875
Concentrate Reconstitution Volume (uL)	875

Calibrator Levels (pg/mL)

A	0.000
B	0.008
C	0.025
D	0.083
E	0.277
F	0.924
G	3.08
H	10.2

Calibration Curve

