

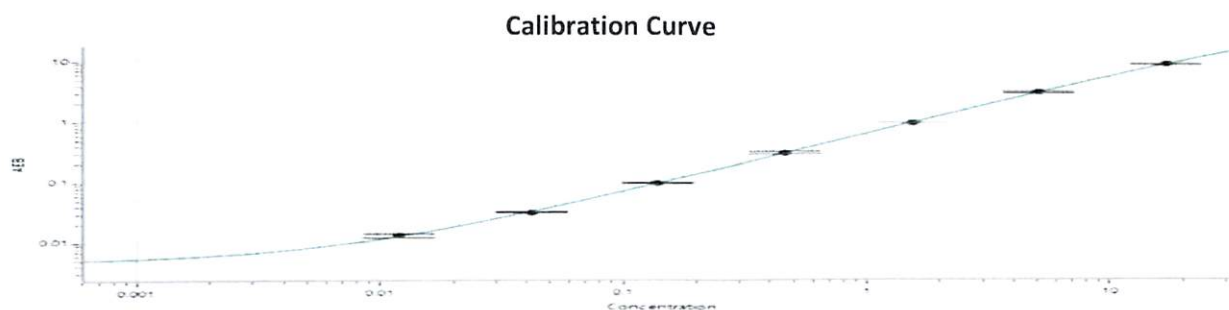
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*

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

**Calibrator Levels
(pg/mL)**

A	0.000
B	0.012
C	0.042
D	0.138
E	0.460
F	1.54
G	5.10
H	17.0



Dhruvi Patel QA Specialist II [Signature] / 09 Dec 2025

Name Title Signature/Date

Release Materials*

IL-13®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.926	0.581 - 1.08
Control 2 Concentrate	412341	22.0	14.8 - 27.4

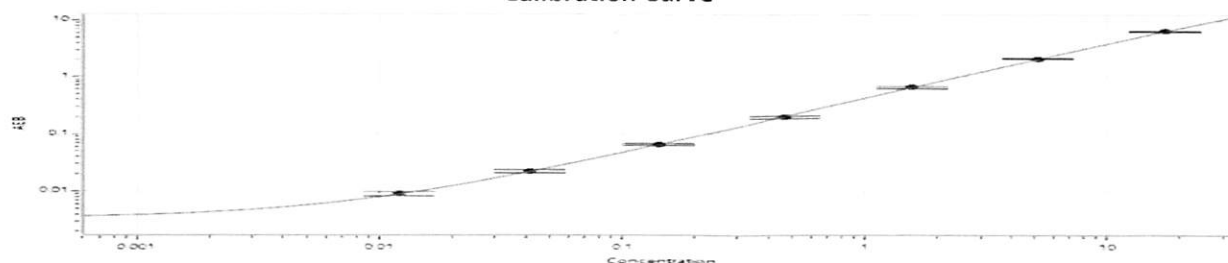
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	0.042
D	0.142
E	0.469
F	1.56
G	5.21
H	17.4

Calibration Curve



Release Materials*

IL-5®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.827	0.549 - 1.02
Control 2 Concentrate	412341	18.3	13.2 - 24.5

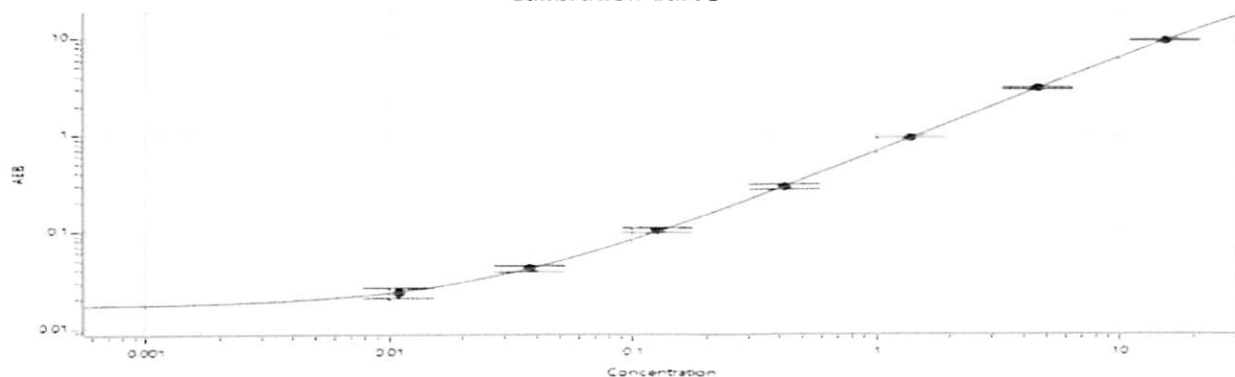
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	0.419
F	1.39
G	4.63
H	15.4

Calibration Curve



Release Materials*

IL-17A®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412340	0.661	0.441 - 0.818
Control 2 Concentrate	412341	14.4	10.3 - 19.1

* 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	0.010
C	0.031
D	0.103
E	0.343
F	1.15
G	3.82
H	12.6

Calibration Curve

