

Product Number:	104979
Lot Number:	504681
Expiration:	29-Jan-2027
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	104964	521910
Detector Reagent	104966	522003
SBG Reagent	104968	522004
Sample Diluent	104969	521012
Calibrator Diluent	104970	521014
Control Diluent	105443	521013
Calibrator Concentrate	104975	411414
Control 1 Concentrate	104977	411416
Control 2 Concentrate	104978	411417
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



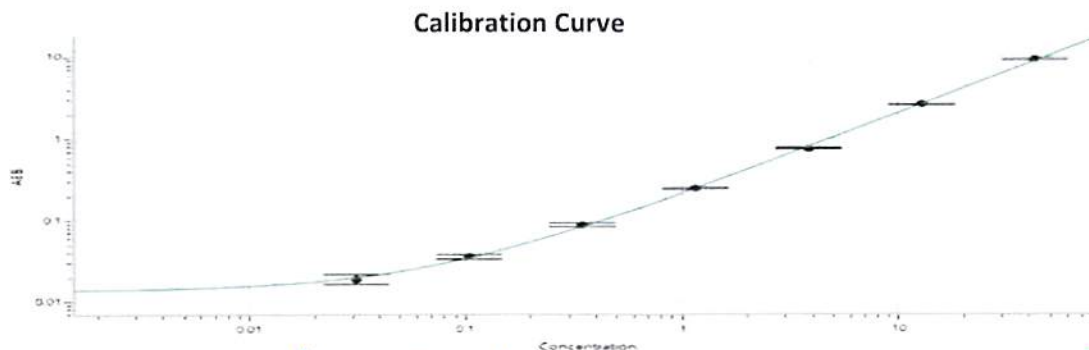
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
IL-6®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	1.97	1.44 - 2.68
Control 2 Concentrate	411417	99.4	70.6 - 131

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

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

Calibrator Levels (pg/mL)	
A	0.000
B	0.031
C	0.104
D	0.345
E	1.15
F	3.83
G	12.7
H	42.5



Review/Approval

Dhruvi Patel
Name

QA Specialist II, QA
Title

SP/24 Sep 2025
Signature/Date

Release Materials*

IL-1Beta®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	0.427	0.252 - 0.588
Control 2 Concentrate	411417	22.6	17.0 - 31.5

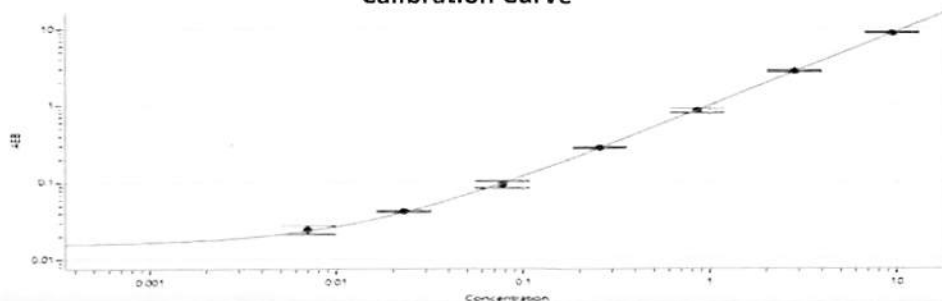
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

Calibrator Levels (pg/mL)

A	0.000
B	0.007
C	0.023
D	0.077
E	0.257
F	0.856
G	2.85
H	9.48

Calibration Curve



Release Materials*

IL-10®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	1.72	0.944 - 2.60
Control 2 Concentrate	411417	76.2	53.9 - 100

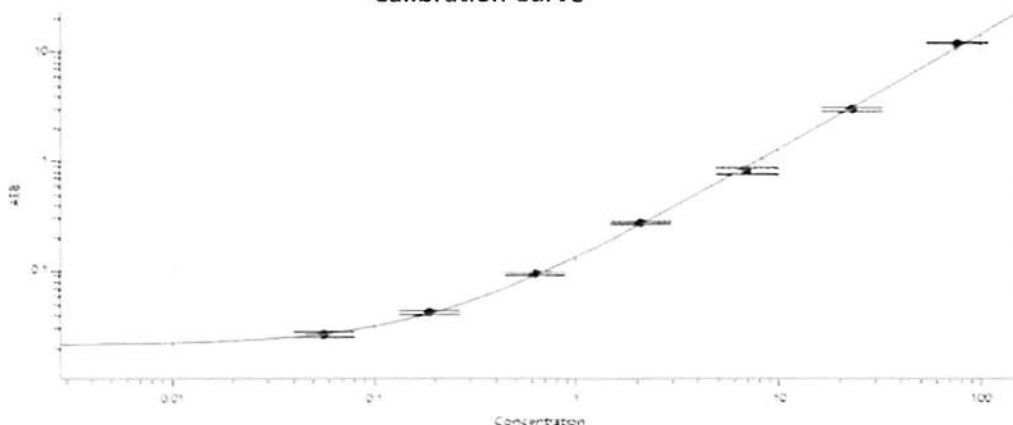
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

Calibrator Levels (pg/mL)

A	0.000
B	0.056
C	0.186
D	0.619
E	2.06
F	6.87
G	22.8
H	76.2

Calibration Curve



Release Materials*

TNFAIpha®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	3.89	2.23 - 5.30
Control 2 Concentrate	411417	175	123 - 230

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

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

**Calibrator Levels
(pg/mL)**

A	0.000
B	0.070
C	0.235
D	0.782
E	2.61
F	8.67
G	28.9
H	96.1

Calibration Curve

