

Product Number:	104979
Lot Number:	504385
Expiration:	7-Feb-2026
Platform(s):	HD-X



Component	Part Number	Lot Number
Bead Reagent	104964	425004
Detector Reagent	104966	425302
SBG Reagent	104968	425303
Sample Diluent	104969	422020
Calibrator Diluent	104970	423321
Control Diluent	105443	423322
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

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

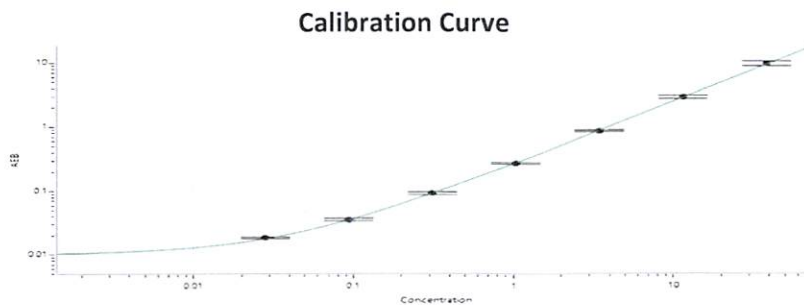
¹ Reagents are not Kit Lot Specific

Release Materials*			
IL-6®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	1.78	1.31 - 2.43
Control 2 Concentrate	411417	91.3	64.6 - 120

* 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	0.028
C	0.094
D	0.312
E	1.04
F	3.46
G	11.5
H	38.4



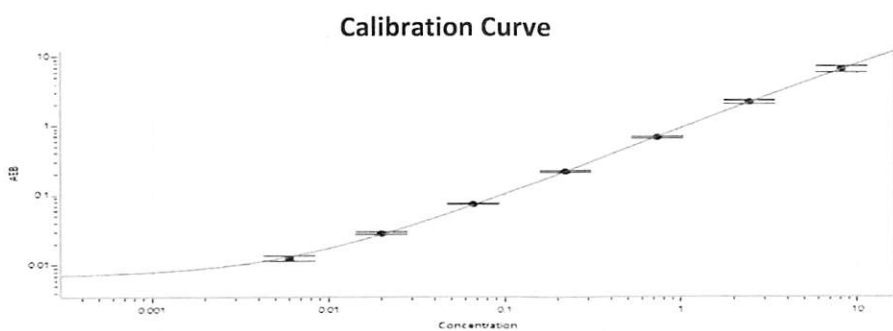
Review/Approval Dhruvi Patel Doc Control I, QA DP 11 Oct 2024
 Name Title Signature/Date

Release Materials*			
IL-1Beta®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	0.388	0.276 - 0.518
Control 2 Concentrate	411417	19.9	14.0 - 26.1

* 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	0.006
C	0.020
D	0.066
E	0.220
F	0.733
G	2.44
H	8.12

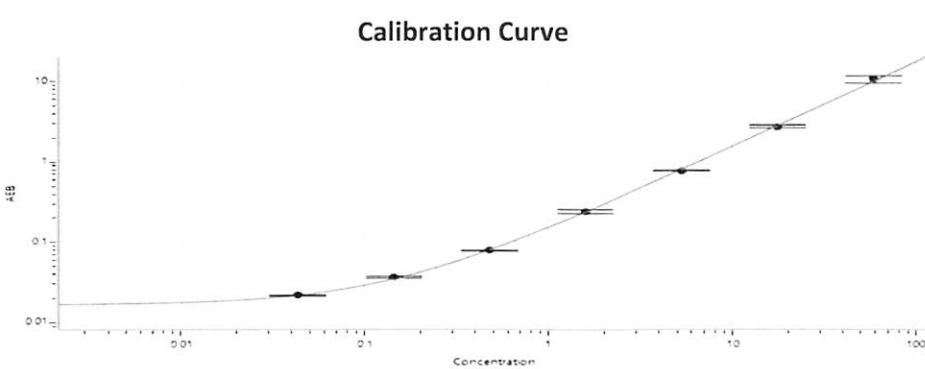


Release Materials*			
IL-10®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	1.29	0.888 - 1.79
Control 2 Concentrate	411417	62.8	45.3 - 84.0

* 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	0.043
C	0.143
D	0.477
E	1.59
F	5.29
G	17.6
H	58.7



Release Materials*

TNFAlpha®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411416	2.84	2.33 - 4.33
Control 2 Concentrate	411417	142	112 - 208

* 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	0.060
C	0.201
D	0.669
E	2.23
F	7.42
G	24.7
H	82.2

Calibration Curve

