

Product Number:	105066
Lot Number:	504419
Expiration:	18-Apr-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	105051	429205	N/A
Detector Reagent	105053	503505	503505A
SBG Reagent	105055	503506	503506A
Sample Diluent	105056	503511	503511A
Calibrator Diluent	105057	430928	N/A
Control Diluent	105398	430927	N/A
Calibrator Concentrate	105062	412220	N/A
Control 1 Concentrate	105064	412221	N/A
Control 2 Concentrate	105065	412222	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

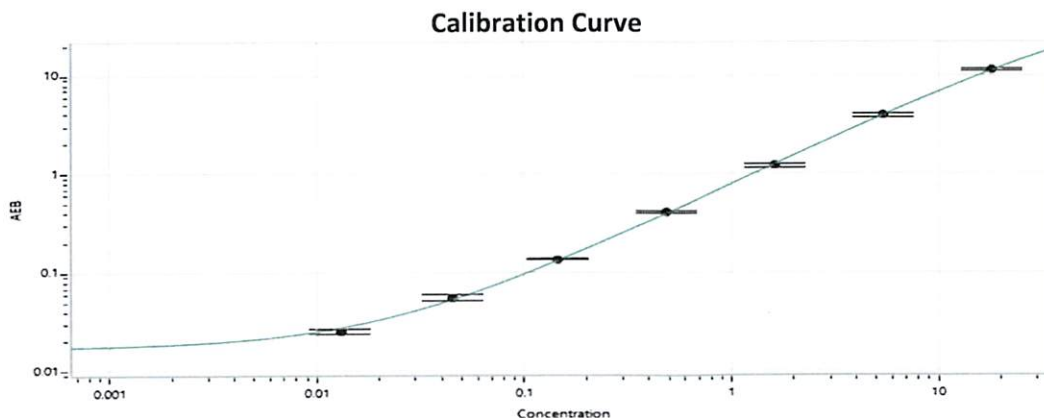
¹ Reagents are not Kit Lot Specific

Release Materials*			
IL-2®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	0.444	0.300 - 0.557
Control 2 Concentrate	412222	20.4	15.1 - 28.0

* 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	0.013
C	0.045
D	0.147
E	0.488
F	1.63
G	5.41
H	18.1



Review/Approval

Dipesh Patel
Name Title

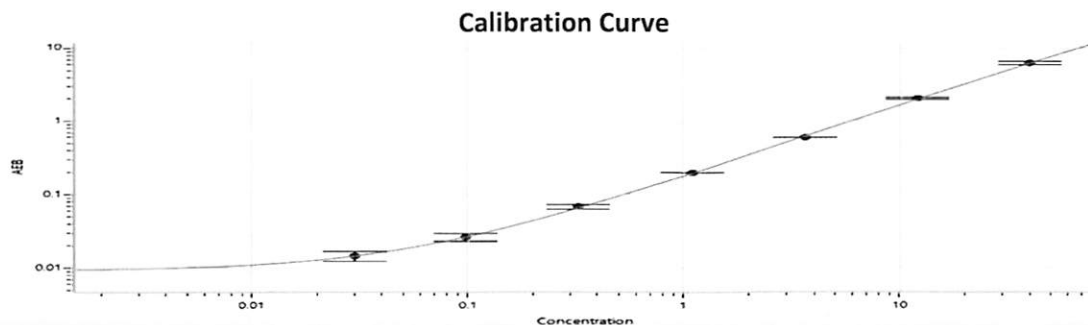
[Signature] / 13 MAR 2025
Signature/Date

Release Materials*			
IFNgamma®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	0.842	0.684 - 1.27
Control 2 Concentrate	412222	45.3	36.9 - 68.5

* 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	0.030
C	0.098
D	0.327
E	1.09
F	3.63
G	12.1
H	40.2

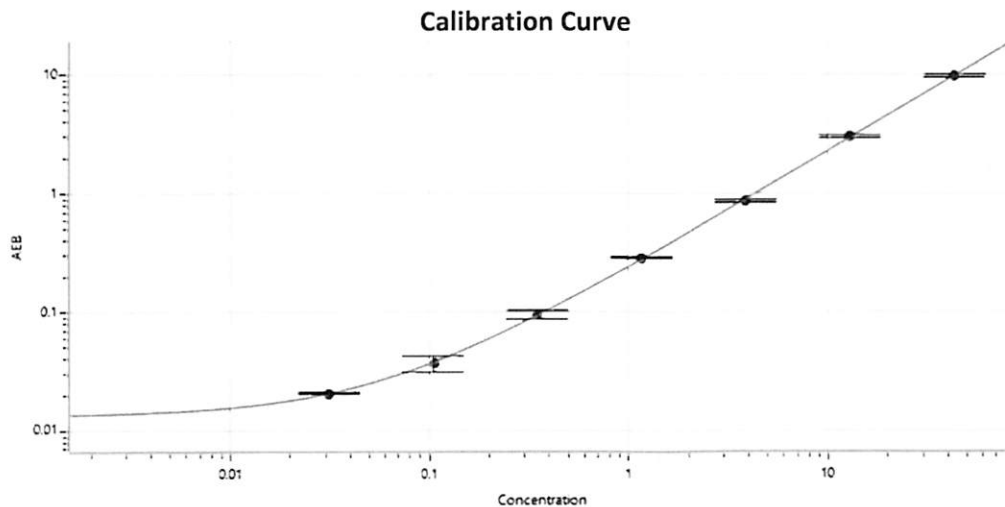


Release Materials*			
IL-6®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	0.925	0.717 - 1.33
Control 2 Concentrate	412222	49.1	34.5 - 64.1

* 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	0.031
C	0.105
D	0.349
E	1.16
F	3.87
G	12.9
H	43.0



Release Materials*

IL-8®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	2.28	1.62 - 3.09
Control 2 Concentrate	412222	114	79.3 - 148

* 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	0.073
C	0.242
D	0.806
E	2.69
F	8.94
G	29.8
H	99.1

Calibration Curve

