

Product Number:	105066
Lot Number:	536616
Expiration:	15-Apr-2027
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	105051	528807
Detector Reagent	105053	529504
SBG Reagent	105055	529505
Sample Diluent	105056	528824
Calibrator Diluent	105057	528825
Control Diluent	105398	528826
Calibrator Concentrate	105062	411437
Control 1 Concentrate	105064	411439
Control 2 Concentrate	105065	411440
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

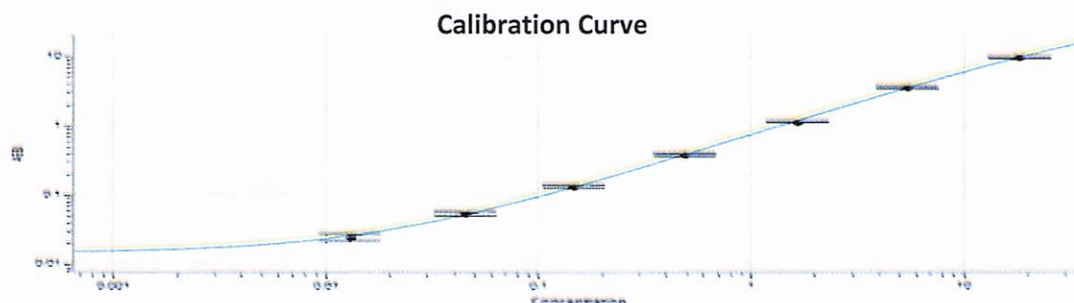
¹ Reagents are not Kit Lot Specific

Release Materials*			
IL-2®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411439	0.489	0.309 - 0.575
Control 2 Concentrate	411440	21.0	15.2 - 28.2

* 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	0.489
F	1.64
G	5.42
H	18.1



Review/Approval

William Chan

Name

Document Control Specialist I

Title

WAC 04DEC2025

Signature/Date

Release Materials*

IFNgamma®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411439	0.892	0.739 - 1.37
Control 2 Concentrate	411440	48.5	34.9 - 64.9

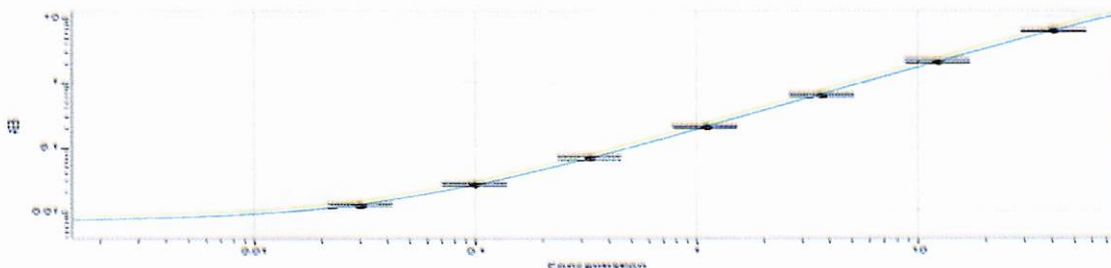
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	3.64
G	12.1
H	40.3

Calibration Curve



Release Materials*

IL-6®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411439	1.17	0.755 - 1.40
Control 2 Concentrate	411440	46.2	34.9 - 64.8

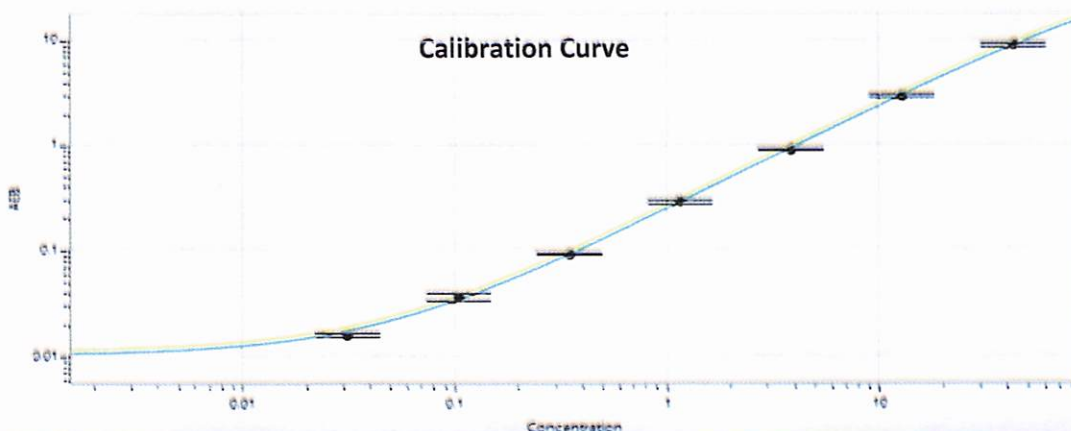
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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

Calibration Curve



IL-8®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	411439	2.81	1.65 - 3.43
Control 2 Concentrate	411440	116	81.6- 152

* 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	0.076
C	0.253
D	0.840
E	2.80
F	9.32
G	31.0
H	103

Calibration Curve

