

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Kit Catalog Numbers: K151QQ-Series
Kit Lot Number: K0081090
Expiration Date: 31 Jan 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-2	80% - 120%	< 20%	90% - 130%*	< 20%	90% - 130%*	< 20%	90% - 130%*	< 20%

* See Additional Comments.

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-2	Calibrator	1420	pg/mL
		Proinflammatory Panel 1 (human) Control 1	390	pg/mL
		Proinflammatory Panel 1 (human) Control 2	38.3	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.34	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1420	543,000	1.4
355	145,000	1.9
88.8	36,900	2.0
22.2	9,010	1.7
5.55	2,270	1.5
1.39	618	1.8
0.347	227	4.2
0.00	93	5.6

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0046914	2-8°C	28 Feb 2021
Anti-hu IL-2 Antibody	D0081129	2-8°C	31 Jan 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080158	2-8°C	31 May 2023
Diluent 2	M0020102	≤ -10°C	28 Feb 2021
Diluent 3	M0030084	≤ -10°C	31 Jan 2021
Read Buffer T	Not Kit Specific	Not Kit Specific	N/A


Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	Not Kit Specific
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory Panel 1 (human) Control 3	Not Kit Specific

Additional Comments:

Assay	Spot	Specification	Comment
1 hu IL-2	1	Control Conc.	Control-1 recovery 130.9%; Control-2 recovery 134.8%, Control-3 recovery 134.8%

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		19 DEC 2018