

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Kit Catalog Numbers: K151QQ-Series; K156QQ-Series;
Kit Lot Number: K0081453
Expiration Date: 31 Jul 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Recovery	CV
		Recovery	CV		
1	IL-2	80% - 120%	< 20%	90% - 130%*	< 20%

* See Additional Comments.

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-2	Calibrator	1420	pg/mL
		Proinflammatory Panel 1 (human) Control 1	384	pg/mL
		Proinflammatory Panel 1 (human) Control 2	37.7	pg/mL
		Proinflammatory Panel 1 (human) Control 3	2.96	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1420	511,000	2.3
355	134,000	5.8
88.8	34,100	1.8
22.2	8,400	4.1
5.55	2,140	1.9
1.39	605	10.0
0.347	202	3.1
0.00	76	5.1

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0047292	2-8°C	31 Aug 2022
Anti-hu IL-2 Antibody	D0081338	2-8°C	31 Jan 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080229	2-8°C	30 Sep 2024
Diluent 2	M0020112	≤ -10°C	31 Jul 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A


Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	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	IL-2	1	Control Conc.	Control-3 recovery 132.2% (specifications 90-130%)

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14 MAY 2020