

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-2 Kit
Kit Catalog Numbers: K152QQ-Series
Kit Lot Number: K0081419
Expiration Date: 31 Mar 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-2	80% - 120%	< 20%	85% - 135%	< 25%	85% - 135%	< 25%	85% - 135%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-2	1.03	75% - 125%	< 25%	1570	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-2	Calibrator	2450	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	919	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	54.8	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	5.84	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
2450	286,000	2.7
613	71,400	1.5
153	18,300	2.4
38.3	4,780	2.4
9.57	1,270	1.9
2.39	359	1.4
0.598	132	3.1
0.00	54	4.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-2 Plate	Z0047261	2-8°C	30 Jun 2022
Anti-ms IL-2 Antibody	D0081345	2-8°C	28 Feb 2023
Proinflammatory panel 1 (mouse) Calibrator Blend	A0080197	2-8°C	31 Dec 2023
Diluent 41	M0410031	≤ -10°C	30 Jun 2022
Diluent 45	M0450031	≤ -10°C	31 Mar 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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

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		23 MAR 2020