

## 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:** K0081537  
**Expiration Date:** 31 Dec 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	2670	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	908	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	54.5	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	5.57	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
2670	382,000	2.3
668	99,300	1.2
167	25,800	2.0
41.7	6,550	1.7
10.4	1,720	2.0
2.61	470	1.4
0.652	165	2.7
0.00	60	4.9

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-2 Plate	Z0047380	2-8°C	28 Feb 2023
Anti-ms IL-2 Antibody	D0081345	2-8°C	28 Feb 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410034	≤ -10°C	31 Dec 2022
Diluent 45	M0450035	≤ -10°C	31 Jan 2023
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
<b>Review/Approval</b>	Karen Hamilla	Quality		05 OCT 2020