

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: K0080950
Expiration Date: 31 Dec 2019

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	ms IL-2	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-2	Calibrator	2870	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	951	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	66.7	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	5.33	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2870	325,000	2.4
718	76,000	2.6
179	17,000	3.6
44.8	4,010	3.6
11.2	1,110	2.9
2.80	337	3.8
0.701	137	4.4
0.00	75	5.1

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-2 Plate	Z0046746	2-8°C	31 Jul 2020
Anti-ms IL-2 Antibody	D0080962	2-8°C	31 Dec 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080138	2-8°C	31 May 2022
Diluent 41	M0410021	≤ -10°C	30 Jun 2020
Diluent 45	M0450020	≤ -10°C	31 May 2020
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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 MAR 2018