



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04440

**Product Description:** Mouse IL-2 Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K152QQD-1, K152QQD-2, K152QQD-4, K152QQG-1, K152QQG-2, K152QQG-4

**Kit Lot Number:** K0080124

**Expiration Date:** 30 JUN 2016

### Acceptance Data:

Metric		Specifications	Assay ms IL-2
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	80% - 130%	Pass
	ULOQ	75% - 125%	Pass
	LLOQ	75% - 125%	Pass
% Recovery CV	Calibrators	< 20%	Pass
	Controls	< 25%	Pass
	ULOQ	< 25%	Pass
	LLOQ	< 25%	Pass

### Concentration of Materials Tested:

Units = pg/mL	Assay ms IL-2
Reconstituted Calibrator	2240
Reconstituted Control 1 (A00C0181)	2630
Reconstituted Control 2 (A00C0182)	193
Reconstituted Control 3 (A00C0183)	17.7
ULOQ	1570
LLOQ	1.03

### Standard Curve Data from Kit QC:

ms IL-2		
Conc., (pg/mL)	Average Signal	%CV
2240	318 078	1.8
560	75 035	3.1
140	19 421	2.6
35.0	4444	5.0
8.75	1264	2.5
2.19	340	2.9
0.547	129	3.5
0	52	4.8



### CERTIFICATE OF ANALYSIS

COA-04440

<b>Mouse IL-2 Kit V-PLEX Components</b>			
Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-2 Plate	Z0045197	2-8°C	31 AUG 2016
Anti-ms IL-2 Antibody	D0080289	2-8°C	31 JUL 2016
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080001	2-8°C	31 DEC 2017
Diluent 41	M0410005	≤ -10°C	30 JUN 2016
Diluent 45	M0450004	≤ -10°C	30 JUN 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
<b>Additional Components Included in V-PLEX Plus</b>			
Wash Buffer (20X), 100 mL	Not Kit Specific	Room Temperature	N/A
Plate Seals	Not Kit Specific	Room Temperature	N/A
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific; See Separate Control Pack COA		
Proinflammatory Panel 1 (mouse) Control 2			
Proinflammatory Panel 1 (mouse) Control 3			

**Statement:** The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dana Short	Quality Control		27 MAY 2014