



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04671

**Product Description:** Mouse IL-6 Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K152QXD-1, K152QXD-2, K152QXD-4, K152QXG-1, K152QXG-2, K152QXG-4

**Kit Lot Number:** K0080202

**Expiration Date:** 30 JUN 2015

### Acceptance Data:

Metric		Specifications	Assay ms IL-6
% 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-6
Reconstituted Calibrator	3710
Reconstituted Control 1 (A00C0181)	5720
Reconstituted Control 2 (A00C0182)	263
Reconstituted Control 3 (A00C0183)	29.2
ULOQ	3140
LLOQ	7.61

### Standard Curve Data from Kit QC:

ms IL-6		
Conc., (pg/mL)	Average Signal	%CV
3710	144 843	4.0
928	28 555	4.2
232	6392	5.4
58.0	1582	3.2
14.5	433	7.8
3.62	165	7.8
0.906	95	7.1
0	78	6.8




## CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-04671

Mouse IL-6 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0044464	2-8°C	30 JUN 2015
Anti-ms IL-6 Antibody	D0080006	2-8°C	30 JUN 2015
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080029	2-8°C	30 SEP 2018
Diluent 41	M0410003	≤ -10°C	31 MAR 2016
Diluent 45	M0450003	≤ -10°C	31 MAR 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
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
Review/Approval	Caitlin Hansen	Quality Control		12 AUG 2014