



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-05300

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: K0080308

Expiration Date: 31 MAR 2017

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	4200
Reconstituted Control 1 (A00C0289)	1740
Reconstituted Control 2 (A00C0290)	271
Reconstituted Control 3 (A00C0291)	66.0
ULOQ	3140
LLOQ	7.61

Standard Curve Data from Kit QC:

ms IL-6		
Conc., (pg/mL)	Average Signal	%CV
4200	125 606	5.0
1050	29 275	6.1
263	6956	5.3
65.6	1759	7.3
16.4	543	12.8
4.10	188	5.1
1.03	102	6.5
0	71	7.1




CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-05300

Mouse IL-6 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0045642	2-8°C	31 AUG 2017
Anti-ms IL-6 Antibody	D0080454	2-8°C	31 MAR 2017
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080043	2-8°C	31 JUL 2019
Diluent 41	M0410008	≤ -10°C	31 JUL 2017
Diluent 45	M0450007	≤ -10°C	31 JUL 2017
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	Dana Short	Quality Control		17 MAR 2015