



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-04428

Product Description: Mouse TNF- α Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K152QWD-1, K152QWD-2, K152QWD-4, K152QWG-1, K152QWG-2, K152QWG-4

Kit Lot Number: K0080145

Expiration Date: 30 APR 2015

Acceptance Data:

Metric		Specifications	Assay ms TNF- α
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	50% - 150%	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 TNF- α
Reconstituted Calibrator	507
Reconstituted Control 1 (A00C0181)	431
Reconstituted Control 2 (A00C0182)	98.0
Reconstituted Control 3 (A00C0183)	21.2
ULOQ	403
LLOQ	0.980

Standard Curve Data from Kit QC:

ms TNF- α		
Conc., (pg/mL)	Average Signal	%CV
507	79 257	4.5
127	15 666	2.9
31.7	4349	2.8
7.92	1077	3.6
1.98	357	3.2
0.495	162	9.7
0.124	115	5.9
0	96	5.7




CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-04428

Mouse TNF-α Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0044472	2-8°C	30 JUN 2015
Anti-ms TNF- α Antibody	D0080020	2-8°C	30 JUN 2016
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080003	2-8°C	31 DEC 2017
Diluent 41	M0410003	$\leq -10^{\circ}\text{C}$	31 MAR 2016
Diluent 45	M0450002	$\leq -10^{\circ}\text{C}$	30 APR 2015
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		12 MAY 2014