



CERTIFICATE OF ANALYSIS

COA-04471

Product Description: Mouse IL-4 Kit V-PLEX® and V-PLEX Plus
Kit Catalog Numbers: K152QRD-1, K152QRD-2, K152QRD-4, K152QRG-1, K152QRG-2, K152QRG-4
Kit Lot Number: K0080140
Expiration Date: 30 APR 2016

Acceptance Data:

Table with 4 columns: Metric, Specifications, Assay, ms IL-4. Rows include % Recovery Concentration and % Recovery CV for Calibrators, Controls, ULOQ, and LLOQ.

Concentration of Materials Tested:

Table with 2 columns: Units = pg/mL, Assay ms IL-4. Rows list Reconstituted Calibrator, Control 1, 2, 3, ULOQ, and LLOQ with their respective values.

Standard Curve Data from Kit QC:

Table with 3 columns: Conc., (pg/mL), Average Signal, %CV. Rows show data points for the standard curve from 1460 down to 0 pg/mL.



CERTIFICATE OF ANALYSIS

COA-04471

<b>Mouse IL-4 Kit V-PLEX Components</b>			
<b>Description</b>	<b>Lot Number</b>	<b>Storage Temperature</b>	<b>Expiration Date</b>
Mouse IL-4 Plate	Z0044454	2-8°C	30 JUN 2015
Anti-ms IL-4 Antibody	D0080014	2-8°C	30 JUN 2015
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080001	2-8°C	31 DEC 2017
Diluent 41	M0410002	≤ -10°C	30 APR 2015
Diluent 45	M0450002	≤ -10°C	30 APR 2015
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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Dana Short	Quality Control		25 MAR 2015




# Meso Scale Diagnostics, L.L.C.

Description: Mouse IL-4 Kit V-PLEX® and V-PLEX Plus

Catalog Number(s): K152QRD/G-1/2/4

Lot #: K0080140

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0080140 has been extended to April 30, 2016.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		24 Mar 2015

This memo supersedes all other statements of date of expiration for this kit, Lot # K0080140.