

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex® and V-Plex Plus Mouse IL-9 Kit
Kit Catalog Numbers: K152XC-Series
Kit Lot Number: K0080745
Expiration Date: 31 Aug 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IL-9	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-9	21.9	75% - 125%	< 20%	2600	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-9	Calibrator	3860	pg/mL
		Cytokine Panel 1 (mouse) Control 1	3670	pg/mL
		Cytokine Panel 1 (mouse) Control 2	893	pg/mL
		Cytokine Panel 1 (mouse) Control 3	209	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-9 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3860	203,000	1.6
965	47,300	3.5
241	10,500	4.0
60.3	2,450	3.0
15.1	692	2.2
3.77	221	2.4
0.942	116	4.6
0.00	74	6.1

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-9 Plate	Z0046391	2-8°C	31 Aug 2019
Anti-ms IL-9 Antibody	D0080934	2-8°C	31 Aug 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080125	2-8°C	31 Mar 2022
Diluent 41	M0410016	2-8°C	31 Aug 2019
Diluent 45	M0450016	2-8°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20x), 100mL	Not Kit Specific
Cytokine Panel 1 (mouse) Control 1	Not Kit Specific
Cytokine Panel 1 (mouse) Control 2	Not Kit Specific
Cytokine Panel 1 (mouse) Control 3	Not Kit Specific

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	23 JUN 2017