

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-5 Kit

Kit Catalog Numbers: K152QS-Series

Kit Lot Number: K0080866

Expiration Date: 31 May 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IL-5	80% - 120%	< 20%	75% - 125%	< 25%	125% - 145%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-5	1.60	75% - 125%	< 25%	590	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-5	Calibrator	969	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	323	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	49.1	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	9.26	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-5 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
969	436,000	1.3
242	103,000	1.5
60.6	23,600	2.4
15.1	6,090	1.6
3.79	1,660	1.7
0.946	480	2.2
0.237	176	1.9
0.00	79	4.8

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-5 Plate	Z0046588	2-8°C	31 Dec 2019
Anti-ms IL-5 Antibody	D0080855	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080138	2-8°C	31 May 2022
Diluent 41	M0410017	≤ -10°C	31 Oct 2019
Diluent 45	M0450018	≤ -10°C	30 Nov 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seal	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory 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		24 OCT 2017