

CERTIFICATE OF ANALYSIS

Page 1 of 2

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

Kit Catalog Numbers: K151QS-Series

Kit Lot Number: K0081001

Expiration Date: 31 Jul 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-5	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-5	6.28	80% - 120%	< 20%	562	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-5	Calibrator	844	pg/mL
		Cytokine Panel 1 (human) Control 1	310	pg/mL
		Cytokine Panel 1 (human) Control 2	110	pg/mL
		Cytokine Panel 1 (human) Control 3	34	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-5 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
844	306,000	1.7
211	78,600	3.0
52.8	19,800	2.8
13.2	5,020	3.5
3.30	1,310	2.7
0.824	374	1.5
0.206	141	3.1
0.00	64	4.9

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-5 Plate	Z0046811	2-8°C	30 Sep 2020
Anti-hu IL-5 Antibody	D0080887	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080121	2-8°C	28 Feb 2022
Diluent 43	M0430024	≤ -10°C	31 Jan 2020
Diluent 3	M0030081	≤ -10°C	30 Sep 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Microplate Adhesive Film	Not Kit Specific
MSD Wash Buffer (20X), 100 mL	Not Kit Specific
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine Panel 1 (human) 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		29 MAY 2018