

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-15 Kit

Kit Catalog Numbers: K151RD-Series

Kit Lot Number: K0081206

Expiration Date: 30 Jun 2021

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-15	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-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-15	Calibrator	829	pg/mL
		Cytokine Panel 1 (human) Control 1	286	pg/mL
		Cytokine Panel 1 (human) Control 2	45	pg/mL
		Cytokine Panel 1 (human) Control 3	8	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
829	284,000	1.4
207	54,800	3.3
51.8	12,700	3.8
13.0	3,100	3.8
3.24	835	2.4
0.810	240	2.1
0.202	110	3.1
0.00	72	7.9

CERTIFICATE OF ANALYSIS

Page 2 of 2

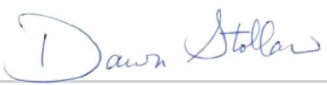
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0047043	2-8°C	31 Aug 2021
Anti-hu IL-15 Antibody	D0081081	2-8°C	30 Sep 2021
Cytokine Panel 1 (human) Calibrator Blend	A0080170	2-8°C	20 Jun 2023
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 2021
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	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		03 APR 2019