

CERTIFICATE OF ANALYSIS

Page 1 of 2

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

Kit Catalog Numbers: K151RD-Series, K156RD-Series

Kit Lot Number: K0081540

Expiration Date: 30 Jun 2022

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-15	Calibrator	941	pg/mL
		Cytokine Panel 1 (human) Control 1	311	pg/mL
		Cytokine Panel 1 (human) Control 2	50.5	pg/mL
		Cytokine Panel 1 (human) Control 3	9.58	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
941	251,000	2.0
235	45,400	2.1
58.8	10,700	2.1
14.7	2,790	2.0
3.68	711	1.9
0.919	213	3.3
0.230	97	5.7
0.00	57	7.6

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0047383	2-8°C	28 Feb 2023
Anti-hu IL-15 Antibody	D0081363	2-8°C	30 Apr 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080240	2-8°C	30 Nov 2024
Diluent 43	M0430037	≤ -10°C	31 Aug 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022
MSD Read Buffer T (4X)	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
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	Karen Hamilla	Quality		09 OCT 2020