

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: K0081036
Expiration Date: 31 Oct 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-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	796	pg/mL
		Cytokine Panel 1 (human) Control 1	288	pg/mL
		Cytokine Panel 1 (human) Control 2	46.3	pg/mL
		Cytokine Panel 1 (human) Control 3	8.87	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
796	234,000	1.8
199	39,200	2.1
49.8	9,350	1.6
12.4	2,350	1.3
3.11	664	1.8
0.777	215	2.9
0.194	99	3.9
0.00	61	5.3

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0046856	2-8°C	31 Dec 2020
Anti-hu IL15 Antibody	D0080983	2-8°C	31 Oct 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080073	2-8°C	31 Jan 2021
Diluent 43	M0430026	≤ -10°C	31 Aug 2020
Diluent 3	M0030082	≤ -10°C	31 Oct 2020
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 (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		22 AUG 2018