

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus® Human IL-7 Kit

Kit Catalog Numbers: K151RC-Series

Kit Lot Number: K0081076

Expiration Date: 30 Apr 2020

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-7	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-7	0.546	80% - 120%	< 20%	563	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-7	Calibrator	844	pg/mL
		Cytokine Panel 1 (human) Control 1	348	pg/mL
		Cytokine Panel 1 (human) Control 2	56.4	pg/mL
		Cytokine Panel 1 (human) Control 3	8.49	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-7 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
844	226,000	3.7
211	54,800	4.9
52.8	12,400	2.4
13.2	3,140	3.4
3.30	800	2.3
0.824	266	4.7
0.206	122	4.7
0.00	76	5.3

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-7 Plate	Z0046901	2-8°C	28 Feb 2021
Anti-hu IL-7 Antibody	D0081055	2-8°C	30 Apr 2020
Cytokine Panel 1 (human) Calibrator Blend	A0080121	2-8°C	28 Feb 2022
Diluent 43	M0430027	≤ -10°C	31 Dec 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 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 (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	<i>Dawn Stollar</i>	04 OCT 2018