

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex® and VPLEX Plus Human IL-3 Kit

Kit Catalog Numbers: K151WZ-Series

Kit Lot Number: K0080677

Expiration Date: 31 Jan 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-3	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-3	12.6	75% - 125%	< 20%	1950	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-3	Calibrator	3360	pg/mL
		Cytokine Panel 2 (human) Control 1	1320	pg/mL
		Cytokine Panel 2 (human) Control 2	282	pg/mL
		Cytokine Panel 2 (human) Control 3	56	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-3 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3360	446,000	4.0
840	64,200	2.3
210	8,960	2.2
52.5	1,660	1.8
13.1	424	2.3
3.28	152	3.1
0.820	84	5.9
0.00	64	5.8

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-3 Plate	Z0046323	2-8°C	31 May 2019
Anti-hu IL-3 Antibody	D0080772	2-8°C	30 Apr 2019
Cytokine Panel 2 (human) Calibrator Blend	A0080114	2-8°C	31 Dec 2021
Diluent 43	M0430018	≤ -10°C	31 Jan 2019
Diluent 3	M0030073	≤ -10°C	31 Mar 2019
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), 100 mL	Not Kit Specific
Cytokine Panel 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (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>	21 JUL 2017