

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 α Kit
Kit Catalog Numbers: K151RB-Series
Kit Lot Number: K0081191
Expiration Date: 28 Feb 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-1 α	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-1 α	2.85	80% - 120%	< 20%	278	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-1 α	Calibrator	454	pg/mL
		Cytokine Panel 1 (human) Control 1	183	pg/mL
		Cytokine Panel 1 (human) Control 2	58.1	pg/mL
		Cytokine Panel 1 (human) Control 3	16.3	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-1 α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
454	71,200	3.1
114	19,300	4.1
28.4	5,120	4.7
7.09	1,350	4.0
1.77	429	3.4
0.443	207	4.5
0.111	154	3.0
0.00	149	13.0

CERTIFICATE OF ANALYSIS

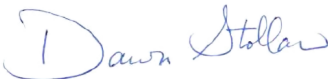
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 α Plate	Z0047028	2-8°C	31 Jul 2021
Anti-hu IL-1 α Antibody	D0081144	2-8°C	28 Feb 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080170	2-8°C	30 Jun 2023
Diluent 43	M0430029	≤ -10°C	30 Apr 2021
Diluent 3	M0030085	≤ -10°C	28 Feb 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), 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	Dawn Stollar	Quality Assurance		11 MAR 2019