

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: K0081585
Expiration Date: 30 Nov 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	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	IL-1 α	2.85	80% - 120%	< 20%	278	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-1 α	Calibrator	454	pg/mL
		Cytokine Panel 1 (human) Control 1	124	pg/mL
		Cytokine Panel 1 (human) Control 2	38.5	pg/mL
		Cytokine Panel 1 (human) Control 3	13	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-1 α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
454	72,500	3.2
114	20,800	3.4
28.4	5,490	2.3
7.09	1,490	2.4
1.77	484	6.9
0.443	239	5.4
0.111	189	19.0
0.00	156	8.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 α Plate	Z0047110	2-8°C	30 Nov 2021
Anti-hu IL-1 α Antibody	D0081144	2-8°C	28 Feb 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080188	2-8°C	30 Nov 2023
Diluent 43	M0430037	≤ -10°C	31 Aug 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022
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	Karen Hamilla	Quality		29 OCT 2020