

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: K0080830
Expiration Date: 31 Jul 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-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	386	pg/mL
		Cytokine Panel 1 (human) Control 1	203	pg/mL
		Cytokine Panel 1 (human) Control 2	61.9	pg/mL
		Cytokine Panel 1 (human) Control 3	20.3	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
386	101,000	1.2
96.5	26,000	2.0
24.1	6,710	1.5
6.03	1,640	1.9
1.51	498	2.6
0.377	178	3.4
0.0942	114	4.8
0.00	149	57.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 α Plate	Z0046543	2-8°C	30 Nov 2019
Anti-hu IL-1 α Antibody	D0080886	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080063	2-8°C	31 Jul 2020
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030077	≤ -10°C	30 Nov 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), 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		02 AUG 2017