

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-1RA Kit
Kit Catalog Numbers: K151WT-Series
Kit Lot Number: K0081224
Expiration Date: 30 Jun 2021

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	IL-1RA	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1RA	9.19	75% - 125%	< 25%	650	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-1RA	Calibrator	732	pg/mL
		Cytokine Panel 2(human) Control 1	514	pg/mL
		Cytokine Panel 2(human) Control 2	93.1	pg/mL
		Cytokine Panel 2(human) Control 3	54.4	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-1RA - Spot 1		
Conc. pg/mL	ECL Signal	%CV
732	221,000	2.1
183	44,600	2.7
45.8	9,570	3.8
11.4	2,420	3.4
2.86	885	18.0
0.715	421	19.0
0.179	397	34.0
0.00	503	36.0

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA Plate	Z0047061	2-8°C	30 Sep 2021
Anti-hu IL-1RA Antibody	D0081093	2-8°C	31 Dec 2021
Cytokine Panel 2(human) Calibrator Blend	A0080126	2-8°C	31 Mar 2022
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 1000mL	Not Kit Specific
Plate Seals	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		09 MAY 2019