

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: K0080689
Expiration Date: 30 Apr 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-1RA	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-1RA	9.19	75% - 125%	< 25%	650	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-1RA	Calibrator	1150	pg/mL
		Cytokine Panel 2 (Human) Control 1	461	pg/mL
		Cytokine Panel 2 (Human) Control 2	81.2	pg/mL
		Cytokine Panel 2 (Human) Control 3	13.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-1RA - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1150	405,000	2.1
288	80,200	2.4
71.9	16,700	3.5
18.0	4,040	7.8
4.49	1,130	5.9
1.12	422	11.0
0.281	237	5.9
0.00	272	16.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1RA Plate	Z0046343	2-8°C	30 Jun 2019
Anti-hu IL-1RA Antibody	D0080771	2-8°C	30 Apr 2019
Cytokine Panel 2 (human) Calibrator Blend	A0080114	2-8°C	31 Dec 2021
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 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), 1000mL	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