

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Rat TNF- α Kit
Kit Catalog Numbers: K153QW-Series
Kit Lot Number: K0080951
Expiration Date: 31 May 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	rt TNF- α	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	rt TNF- α	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	rt TNF- α	Calibrator	1360	pg/mL
		Proinflammatory Panel 1 (rat) Control 1	1200	pg/mL
		Proinflammatory Panel 1 (rat) Control 2	357	pg/mL
		Proinflammatory Panel 1 (rat) Control 3	102	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

rt TNF- α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1360	310,000	1.4
340	59,700	0.9
85.0	13,000	1.6
21.3	3,100	1.7
5.31	874	2.2
1.33	292	3.4
0.332	139	4.2
0.00	90	5.9

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat TNF- α Plate	Z0046747	2-8°C	31 Jul 2020
Anti-rat TNF- α Antibody	D0081033	2-8°C	29 Feb 2020
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080124	2-8°C	31 Mar 2022
Blocker H	R0100015	Room Temperature	30 Sep 2019
Diluent 42	M0420010	$\leq -10^{\circ}\text{C}$	31 Jul 2020
Diluent 40	M0400010	$\leq -10^{\circ}\text{C}$	31 May 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
Proinflammatory Panel 1 (rat) Control 1	Not Kit Specific
Proinflammatory Panel 1 (rat) Control 2	Not Kit Specific
Proinflammatory Panel 1 (rat) 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>	17 APR 2018