

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: K0081417
Expiration Date: 31 Aug 2021

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	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	TNF- α	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	TNF- α	Calibrator	1200	pg/mL
		Proinflammatory Panel 1 (rat) Control 1	948	pg/mL
		Proinflammatory Panel 1 (rat) Control 2	293	pg/mL
		Proinflammatory Panel 1 (rat) Control 3	113	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

TNF- α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1200	238,000	4.3
300	47,300	4.8
75.0	9,960	4.5
18.8	2,400	5.5
4.69	718	5.0
1.17	241	3.6
0.293	130	6.6
0.00	90	6.7

Components:

CERTIFICATE OF ANALYSIS


Page 2 of 2

Description	Lot Number	Storage Temperature	Expiration Date
Rat TNF- α Plate	Z0047267	2-8°C	31 Jul 2022
Anti-rat TNF- α Antibody	D0081204	2-8°C	31 May 2022
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080177	2-8°C	31 Dec 2023
Blocker H	R0100018	Room Temperature	31 Mar 2022
Diluent 42	M0420012	$\leq -10^{\circ}\text{C}$	30 Sep 2021
Diluent 40	M0400014	$\leq -10^{\circ}\text{C}$	31 Aug 2021
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
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	Karen Hamilla	Quality		27 MAR 2020