

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse TNF- α Kit
Kit Catalog Numbers: K152QW- Series
Kit Lot Number: K0081319
Expiration Date: 31 Dec 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) 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%

* See Additional Comments.

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- α	0.980	75% - 125%	< 25%	403	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	TNF- α	Calibrator	629	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	166	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	29.9	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.93	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
629	120,000	2.2
157	27,600	3.1
39.3	6,890	2.2
9.83	1,730	2.0
2.46	476	2.6
0.614	180	2.9
0.154	108	3.6
0.00	82	5.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0047167	2-8°C	28 Feb 2022
Anti-ms TNF-a Antibody	D0081226	2-8°C	31 May 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080168	2-8°C	31 May 2023
Diluent 41	M0410029	$\leq -10^{\circ}\text{C}$	31 Dec 2021
Diluent 45	M0450030	$\leq -10^{\circ}\text{C}$	28 Feb 2022


Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X)	Not Kit Specific
Plate Seals	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

Additional Comments:

	Assay	Spot	Specification	Comment
1	TNF- α	1	Control % Recovery	Control 1 = 136.2% Control 2 = 130.5%

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		18 NOV 2019