

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: K0081539

Expiration Date: 31 May 2022

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%

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	516	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	158	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	28.1	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.46	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
516	125,000	2.2
129	28,500	2.0
32.3	6,960	2.2
8.06	1,730	1.7
2.02	486	3.3
0.504	186	3.8
0.126	116	4.6
0.00	91	4.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0047382	2-8°C	28 Feb 2023
Anti-ms TNF- α Antibody	D0081226	2-8°C	31 May 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410034	\leq -10°C	31 Dec 2022
Diluent 45	M0450035	\leq -10°C	31 Jan 2023
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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) 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		08 OCT 2020