

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: K0081075
Expiration Date: 30 Jun 2020

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	ms 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	ms TNF- α	0.980	75% - 125%	< 25%	403	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms TNF- α	Calibrator	662	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:

ms TNF- α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
662	159,000	2.6
166	35,700	2.5
41.4	8,310	1.6
10.3	2,030	1.4
2.59	577	2.8
0.646	213	5.6
0.162	123	11.0
0.00	96	9.2

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0046900	2-8°C	28 Feb 2021
Anti-ms TNF- α Antibody	D0081060	2-8°C	30 Jun 2020
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080152	2-8°C	30 Nov 2022
Diluent 41	M0410024	\leq -10°C	30 Sep 2020
Diluent 45	M0450021	\leq -10°C	30 Sep 2020
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

Additional Comments:

Assay	Spot	Specification	Comment
1 ms TNF- α	1	Control Conc.	Control-1 recovery 136.6%; Control-2 recovery 130%

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		08 OCT 2018