

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: K0081153
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	145,000	3.3
166	34,200	5.7
41.4	7,790	5.0
10.3	2,070	4.7
2.59	568	4.0
0.646	215	3.8
0.162	113	4.4
0.00	91	8.2

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0046976	2-8°C	30 Apr 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	M0410025	$\leq -10^{\circ}\text{C}$	31 Jan 2021
Diluent 45	M0450023	$\leq -10^{\circ}\text{C}$	31 Jan 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 (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-2 recovery : 125.3%

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>	14 DEC 2018