

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: K0081079
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	Value	LLOQ			Value	ULOQ	
			Recovery	CV	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	674	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
674	162,000	2.4
169	36,400	2.6
42.1	8,880	2.0
10.5	2,260	4.0
2.63	607	2.8
0.658	214	4.3
0.165	122	5.6
0.00	94	5.4

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0046764	2-8°C	31 Jul 2020
Anti-ms TNF- α Antibody	D0081060	2-8°C	30 Jun 2020
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080138	2-8°C	31 May 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

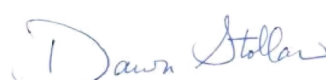
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

Additional Comments:

Assay	Spot	Specification	Comment
1 ms TNF- α	1	Control Conc.	% Recovery for Control 1 = 128.3%

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		15 APR 2019