

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: K0080836
Expiration Date: 31 Mar 2019

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	653	pg/mL
		Proinflammatory Panel 1 (Mouse) Control 1	140	pg/mL
		Proinflammatory Panel 1 (Mouse) Control 2	27.8	pg/mL
		Proinflammatory Panel 1 (Mouse) Control 3	5.12	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
653	194,000	2.7
163	45,700	1.8
40.8	10,900	2.9
10.2	2,780	2.5
2.55	815	2.0
0.638	288	3.5
0.159	161	2.8
0.00	116	3.4

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0046554	2-8°C	30 Nov 2019
Anti-ms TNF- α Antibody	D0080787	2-8°C	31 Mar 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080110	2-8°C	31 Dec 2021
Diluent 41	M0410017	\leq -10°C	31 Oct 2019
Diluent 45	M0450018	\leq -10°C	30 Nov 2019
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.	The observed % Recovery for Control 1 was 139.8% and for Control 2 was 135.3%.

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	31JUL2017