



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse TNF- α Kit
Kit Lot Number: K0081982
Expiration Date: 30 Sep 2023
Kit Catalog Numbers: K152QW-Series

Precision and Accuracy 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%

* See Additional Comments.

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 Calibrator:

Spot	Name	Conc.	Units
1	TNF- α	516	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse TNF- α Plate	Z0047670	2-8°C	31 Aug 2024
Anti-ms TNF- α Antibody	D0081489	2-8°C	30 Sep 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410043	\leq -10°C	31 Mar 2024
Diluent 45	M0450044	\leq -10°C	31 Mar 2024

Additional Comments:

Assay	Spot	Specification	Comment
TNF- α	1	Control Conc.	% Control recovery specification 125-150%, the actual measured % recovery was 133.8%

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

Name	Function	Signature	Date
Review/Approval	Quality		18 JUL 2022