

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Rat TNF- $\alpha$  Kit  
**Kit Lot Number:** K0081509  
**Expiration Date:** 31 Aug 2022  
**Kit Catalog Numbers:** K153QW-Series, K153B2-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- $\alpha$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF- $\alpha$	9.10	75% - 125%	< 25%	793	75% - 125%	< 25%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	TNF- $\alpha$	1200	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat TNF- $\alpha$ Plate	Z0047355	2-8°C	31 Dec 2022
Anti-rat TNF- $\alpha$ Antibody	D0081472	2-8°C	31 Aug 2023
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080177	2-8°C	31 Dec 2023
Blocker H	R0100019	Room Temperature	31 Aug 2022
Diluent 42	M0420013	$\leq -10^{\circ}\text{C}$	31 Dec 2022
Diluent 40	M0400016	$\leq -10^{\circ}\text{C}$	30 Nov 2022

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	Karen Hamilla	Quality		01 FEB 2021