

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Rat TNF- α Kit
Kit Lot Number: K0081957
Expiration Date: 30 Apr 2023
Kit Catalog Numbers: K153QW-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- α	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	TNF- α	1230	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat TNF- α Plate	Z0047645	2-8°C	30 Jun 2024
Anti-rat TNF- α Antibody	D0081641	2-8°C	31 Aug 2024
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080290	2-8°C	30 Apr 2026
Blocker H	R0100020	2-8°C	30 Apr 2023
Diluent 42	M0420016	\leq -10°C	31 Jul 2024
Diluent 40	M0400018	\leq -10°C	31 Jul 2024

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		13 JUL 2022