

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human TNF- α Kit
Kit Lot Number: K0081931
Expiration Date: 30 Sep 2023
Kit Catalog Numbers: K151QW-Series

Precision and Accuracy Specifications:

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TNF- α	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	TNF- α	0.69	80% - 120%	< 20%	248	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	TNF- α	348	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- α Plate	Z0047638	2-8°C	30 Jun 2024
Anti-hu TNF- α Antibody	D0081777	2-8°C	31 Aug 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080310	2-8°C	31 Oct 2026
Diluent 2	M0020116	\leq -10°C	30 Sep 2023
Diluent 3	M0030111	\leq -10°C	31 May 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		31 MAR 2022