

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-Plex Plus Human TNF- α Kit
Kit Catalog Numbers: K151QW-Series
Kit Lot Number: K0081310
Expiration Date: 31 Dec 2021

Acceptance Data/ Percent Recovery Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TNF-a	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	TNF-a	Calibrator	320	pg/ml
		Proinflammatory Panel 1 (human) Control 1	87.8	pg/ml
		Proinflammatory Panel 1 (human) Control 2	15.7	pg/ml
		Proinflammatory Panel 1 (human) Control 3	2.02	pg/ml

Controls tested may not be the same lot as shipped.

Standard Curve Data:

TNF-a - Spot 1		
Conc. pg/ml	ECL Signal	%CV
320	124,000	2.1
80.0	30,300	2.5
20.0	7,770	2.8
5.00	1,980	1.9
1.25	546	2.0
0.313	183	2.8
0.0781	92	4.4
0.00	59	7.5

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TNF- α Plate	Z0047155	2-8°C	28 Feb 2022
Anti-hu TNF- α Antibody	D0081308	2-8°C	31 Oct 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080216	2-8°C	20 Jun 2024
Diluent 2	M0020109	\leq -10°C	28 Feb 2022
Diluent 3	M0030090	\leq -10°C	31 Dec 2021
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 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory Panel 1 (human) Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 DEC 2019