

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Proinflammatory Panel1 (human) Kit**Lot Number:** K0082993**Expiration Date:** 31 JUL 2028**Catalog Number:** K15049-Series, K15052-Series, K15053-Series, K15056-Series, K151QR-Series, K151OD-Series, K15345-Series, K15348-Series, K15346-Series, K15349-Series, K156AR-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- γ Antibody	D0082629	2-8°C	31 OCT 2029
Anti-hu IL-10 Antibody	D0082505	2-8°C	31 JAN 2029
Anti-hu IL-12p70 Antibody	D0082433	2-8°C	31 AUG 2028
Anti-hu IL-13 Antibody	D0082514	2-8°C	28 FEB 2029
Anti-hu IL-1 β Antibody	D0082597	2-8°C	30 SEP 2029
Anti-hu IL-2 Antibody	D0082621	2-8°C	31 AUG 2029
Anti-hu IL-4 Antibody	D0082598	2-8°C	31 JUL 2029
Anti-hu IL-6 Antibody	D0082569	2-8°C	30 JUN 2029
Anti-hu IL-8 Antibody	D0082504	2-8°C	31 JAN 2029
Anti-hu TNF- α Antibody	D0082624	2-8°C	31 AUG 2029
Proinflammatory Panel 1 (hu) 96-Well 10-Spot SECTOR Plate	Z0048283	2-8°C	31 DEC 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080464	2-8°C	30 APR 2031
Diluent 2	M0020138	$\leq -10^{\circ}\text{C}$	30 SEP 2028
Diluent 3	M0030148	$\leq -10^{\circ}\text{C}$	31 JUL 2028

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

Precision, Accuracy, and Sensitivity 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	IFN- γ	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	IL-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	IL-2	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	IL-4	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
5	IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
6	IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	IL-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	IL-12p70	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
9	IL-13	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	TNF- α	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%



CERTIFICATE OF ANALYSIS

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%
2	IL-1 β	0.646	80% - 120%	< 20%	375	80% - 120%	< 20%
3	IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%
4	IL-4	0.218	80% - 120%	< 20%	158	80% - 120%	< 20%
5	IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%
6	IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%
7	IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%
8	IL-12p70	1.22	80% - 120%	< 20%	315	80% - 120%	< 20%
9	IL-13	4.21	80% - 120%	< 20%	353	80% - 120%	< 20%
10	TNF- α	0.690	80% - 120%	< 20%	248	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	Spec
1	IFN- γ	< 1.0%
2	IL-1 β	< 1.0%
3	IL-2	< 1.0%
4	IL-4	< 1.0%
5	IL-6	< 1.0%
6	IL-8	< 1.0%
7	IL-10	< 1.0%
8	IL-12p70	< 1.0%
9	IL-13	< 1.0%
10	TNF- α	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	1520	pg/mL
2	IL-1 β	595	pg/mL
3	IL-2	1720	pg/mL
4	IL-4	254	pg/mL
5	IL-6	796	pg/mL
6	IL-8	583	pg/mL
7	IL-10	375	pg/mL
8	IL-12p70	479	pg/mL
9	IL-13	531	pg/mL
10	TNF- α	344	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 AUG 2026