



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Proinflammatory Panel1 (human) Kit

Lot Number: K0082924

Expiration Date: 31 DEC 2027

Catalog Number: K15049-Series, K15052-Series, K15053-Series, K15056-Series, K151QR-Series, K151OD-Series, K15345-Series, K15348-Series, K15346-Series, K15349-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	D0082502	2-8°C	31 JAN 2029
Anti-hu IL-10 Antibody	D0082432	2-8°C	31 AUG 2028
Anti-hu IL-12p70 Antibody	D0082529	2-8°C	31 MAR 2029
Anti-hu IL-13 Antibody	D0082471	2-8°C	31 OCT 2028
Anti-hu IL-1 β Antibody	D0082597	2-8°C	30 SEP 2029
Anti-hu IL-2 Antibody	D0082470	2-8°C	31 OCT 2028
Anti-hu IL-4 Antibody	D0082287	2-8°C	31 DEC 2027
Anti-hu IL-6 Antibody	D0082503	2-8°C	31 JAN 2029
Anti-hu IL-8 Antibody	D0082431	2-8°C	31 AUG 2028
Anti-hu TNF- α Antibody	D0082515	2-8°C	28 FEB 2029
Proinflammatory Panel 1 (hu) 96-Well 10-Spot SECTOR Plate	Z0048248	2-8°C	30 SEP 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080442	2-8°C	31 AUG 2030
Diluent 2	M0020136	$\leq -10^{\circ}\text{C}$	31 JAN 2028
Diluent 3	M0030147	$\leq -10^{\circ}\text{C}$	31 MAY 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

Spot	Name	LLOQ			ULOQ		
		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	1600	pg/mL
2	IL-1 β	562	pg/mL
3	IL-2	1510	pg/mL
4	IL-4	245	pg/mL
5	IL-6	730	pg/mL
6	IL-8	557	pg/mL
7	IL-10	348	pg/mL
8	IL-12p70	457	pg/mL
9	IL-13	517	pg/mL
10	TNF- α	346	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		29 APR 2026