

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX and V-PLEX Plus Proinflammatory Panel1 (human) Kit
Lot Number: K0082898
Expiration Date: 30 SEP 2027
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	D0082502	2-8°C	31 JAN 2029
Anti-hu IL-10 Antibody	D0082251	2-8°C	30 SEP 2027
Anti-hu IL-12p70 Antibody	D0082433	2-8°C	31 AUG 2028
Anti-hu IL-13 Antibody	D0082471	2-8°C	31 OCT 2028
Anti-hu IL-1 β Antibody	D0082512	2-8°C	30 JUN 2029
Anti-hu IL-2 Antibody	D0082470	2-8°C	31 OCT 2028
Anti-hu IL-4 Antibody	D0082441	2-8°C	31 AUG 2028
Anti-hu IL-6 Antibody	D0082503	2-8°C	31 JAN 2029
Anti-hu IL-8 Antibody	D0082504	2-8°C	31 JAN 2029
Anti-hu TNF- α Antibody	D0082515	2-8°C	28 FEB 2029
Proinflammatory Panel 1 (hu) 96-Well 10-Spot SECTOR Plate	Z0048236	2-8°C	31 AUG 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080439	2-8°C	31 JUL 2030
Diluent 2	M0020136	$\leq -10^{\circ}\text{C}$	31 JAN 2028
Diluent 3	M0030144	$\leq -10^{\circ}\text{C}$	31 JAN 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	0.89	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	0.69	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	1815	pg/mL
2	IL-1 β	592	pg/mL
3	IL-2	1580	pg/mL
4	IL-4	260	pg/mL
5	IL-6	783	pg/mL
6	IL-8	562	pg/mL
7	IL-10	366	pg/mL
8	IL-12p70	486	pg/mL
9	IL-13	547	pg/mL
10	TNF- α	365	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	Michelle Likiam	Quality		24 MAR 2026