



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human TARC Gen. B Kit, SECTOR
Lot Number: K0082933
Expiration Date: 31 MAY 2027
Catalog Number: K151ANK-Series, K156ANK-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
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080416	2-8°C	31 DEC 2029
Human TARC Antibody, 1 Plate	D0082398	2-8°C	30 JUN 2028
Human TARC Gen. B 96-Well Small Spot SECTOR Plate	Z0048253	2-8°C	30 SEP 2028
Diluent 3	M0030137	≤ -10°C	31 JUL 2027
Diluent 57	M0570018	≤ -10°C	31 MAY 2027

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Gen. B Control 1		Chemokine Panel 1 (human) Gen. B Control 2		Chemokine Panel 1 (human) Gen. B Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TARC/CCL17	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TARC/CCL17	2.14	80% - 120%	< 20%	481	75% - 125%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TARC/CCL17	744	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		13 MAY 2026