



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IL-15 Kit**Lot Number:** K0082955**Expiration Date:** 28 FEB 2027**Catalog Number:** K151RD-Series, K156RD-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
Human IL-15, 96-Well Small Spot SECTOR Plate	Z0048171	2-8°C	29 FEB 2028
Cytokine Panel 1 (human) Calibrator Blend	A0080454	2-8°C	31 DEC 2030
Anti-hu IL-15 Antibody	D0082181	2-8°C	30 JUN 2027
Diluent 3	M0030140	≤ -10°C	30 SEP 2027
Diluent 43	M0430058	≤ -10°C	28 FEB 2027

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-15	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-15	846	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		08 JUL 2026