



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IFN- γ V2 Kit, SECTOR

Lot Number: K0082728

Expiration Date: 31 JUL 2027

Catalog Number: K151AVC-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Antibody	D0082498	2-8°C	31 DEC 2028
Human IFN- γ V2 96-Well Small Spot SECTOR Plate	Z0048135	2-8°C	30 NOV 2027
Diluent 58	M0580015	\leq -10°C	31 JUL 2027
Cytolytic Panel 1 (human) Calibrator Blend	A0080420	\leq -70°C	31 JUL 2027

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Cytolytic Panel 1 (human) Control 1		Cytolytic Panel 1 (human) Control 2		Cytolytic 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%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.748	80% - 120%	< 20%	1100	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	29700	pg/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		16 SEP 2025