

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-Plex Plus Human IL-2 Kit
Kit Lot Number: K0081627
Expiration Date: 31 May 2023
Kit Catalog Numbers: K151QQ-Series, K156QQ-Series

Precision and Accuracy 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	IL-2	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-2	1500	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0047450	2-8°C	30 Jun 2023
Anti-hu IL-2 Antibody	D0081510	2-8°C	30 Sep 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080281	2-8°C	28 Feb 2026
Diluent 2	M0020115	≤ -10°C	31 May 2023
Diluent 3	M0030106	≤ -10°C	31 Aug 2023

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 APR 2021