

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-Plex® and V-Plex Plus Mouse IL-2 Kit
Kit Lot Number: K0081675
Expiration Date: 28 Feb 2023
Kit Catalog Numbers: K152QQ-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-2	80% - 120%	< 20%	85% - 135%	< 25%	85% - 135%	< 25%	85% - 135%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-2	1.03	75% - 125%	< 25%	1570	75% - 125%	< 25%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-2	2670	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-2 Plate	Z0047475	2-8°C	31 Jul 2023
Anti-ms IL-2 Antibody	D0081345	2-8°C	28 Feb 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410039	≤ -10°C	31 Aug 2023
Diluent 45	M0450039	≤ -10°C	30 Jun 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		26 MAY 2021