



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Mouse IL-2 Kit


Catalog Number(s) / Size(s): K152QQ - Series

Lot Number: K0081810

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081810 has been extended to 29 FEB 2024.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse IL-2 Plate	Z0047475	31 JUL 2023	29 FEB 2024
Anti-ms IL-2 Antibody	D0081345	28 FEB 2023	29 FEB 2024
Diluent 41	M0410041	30 NOV 2023	29 FEB 2024
Diluent 45	M0450042	31 DEC 2023	29 FEB 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		11 JAN 2023

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-2 Kit
Kit Lot Number: K0081810
Expiration Date: 28 Feb 2023
Kit Catalog Numbers: K152QQ-Series

Precision and Accuracy Specifications:

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

		LLOQ			ULOQ		
Spot	Name	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	M0410041	≤ -10°C	30 Nov 2023
Diluent 45	M0450042	≤ -10°C	31 Dec 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		09 AUG 2021