



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Mouse MIP-3 α Kit

Catalog Number(s) / Size(s): K152XD - Series

Lot Number: K0080775

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080775 has been extended to 30 SEP 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse MIP-3 α Plate	Z0046459	30 SEP 2019	30 SEP 2020
Anti-ms MIP-3 α Antibody	D0080949	31 OCT 2019	30 SEP 2020
Diluent 41	M0410018	30 NOV 2019	30 SEP 2020
Diluent 45	M0450019	31 DEC 2019	30 SEP 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		28 FEB 2020

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

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse MIP-3 α Kit
Kit Catalog Numbers: K152XD-Series
Kit Lot Number: K0080775
Expiration Date: 30 Sep 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (mouse) Control 1		TH17 Panel 1 (mouse) Control 2		TH17 Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms MIP-3 α	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms MIP-3 α	3.42	75% - 125%	< 20%	530	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms MIP-3 α	Calibrator	1130	pg/mL
		TH17 Panel 1 (mouse) Control 1	788	pg/mL
		TH17 Panel 1 (mouse) Control 2	188	pg/mL
		TH17 Panel 1 (mouse) Control 3	56.1	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms MIP-3 α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1130	344,000	2.0
283	66,900	3.5
70.6	13,700	3.9
17.7	3,530	3.2
4.41	905	3.2
1.10	285	6.1
0.276	124	3.1
0.00	71	6.9

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse MIP-3 α Plate	Z0046459	2-8°C	30 Sep 2019
Anti-ms MIP-3 α Antibody	D0080949	2-8°C	31 Oct 2019
TH17 Panel 1 (mouse) Calibrator Blend	A0080146	2-8°C	31 Aug 2022
Diluent 41	M0410018	\leq -10°C	30 Nov 2019
Diluent 45	M0450019	\leq -10°C	31 Dec 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100 mL	Not Kit Specific
TH17 Panel 1 (mouse) Control 1	Not Kit Specific
TH17 Panel 1 (mouse) Control 2	Not Kit Specific
TH17 Panel 1 (mouse) Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	02 JAN 2018