



RFE MEMO

Page 1 of 1

Product Description(s): MSD® Mouse/Rat Insulin Kit

Catalog Number(s) / Size(s): K152BZC-Series

Lot Number: K0014046

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0014046 has been extended to 30 Apr 2020.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse/Rat Insulin Plate	Z0055932	31 OCT 2019	30 APR 2020
Insulin Calibrator	A0010403	30 APR 2019	30 APR 2020
Anti-m/r insulin Antibody	D0013654	31 JUL 2019	30 APR 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		03 SEP 2019

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



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Mouse/Rat Insulin Kit

Catalog Number: K152BZC - Series

Lot Number: K0014046

Expiration Date: N/A (See expiration date on specific components below)

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Kit Contents:

Plate					
Lot #	Z0055932	Catalog #	L452BZA	Expiration Date:	31 OCT 2019

Antibody Blend	
Lot Number:	D0013654
Catalog Number:	D22BZ
Expiration Date:	31 JUL 2019
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010403
Catalog Number:	C00BZ
Expiration Date:	30 APR 2019
Storage Temperature:	≤ -70°C

Specifications:

Release method:

Coating Uniformity	Specifications
Mean Intraplate CV	≤ 15%
Mean Interplate CV	≤ 18%
No more than one plate with CV	> 18%
No plate with CV	> 20%

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>	08 NOV 2017