



RFE MEMO

Page 1 of 1

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

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

Lot Number: K0014172

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0014172 has been extended to 31 JAN 2021.

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	Z0056028	31 JAN 2020	31 JAN 2021
Anti-m/r Insulin Antibody	D0013688	31 JAN 2020	31 JAN 2021
Insulin Calibrator	A0010433	31 APR 2020	31 JAN 2021

Function	Name	Signature	Date
Quality	Karen Hamilla		04 DEC 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: K0014172

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 #	Z0056028	Catalog #	L452BZA	Expiration Date:	31 JAN 2020

Antibody Blend	
Lot Number:	D0013688
Catalog Number:	D22BZ
Expiration Date:	31 JAN 2020
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0010433
Catalog Number:	C00BZ
Expiration Date:	30 APR 2020
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>	02 APR 2018