



CERTIFICATE OF ANALYSIS

COA - 07854

Product Description: MSD[®] Mouse Metabolic Kit

Catalog #/Size: K15124C - Series

Lot Number: K0013971

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 Number: Z0055682

Plate Lot Catalog Number: N45124A

Plate Lot Expiration Date: 31 JAN 2019

Antibody blend Lot Number: D0013582

Antibody blend Catalog Number: D2124

Antibody blend Expiration Date: 31 JAN 2019

Calibrator blend Lot Number: A0010397

Calibrator blend Catalog Number: C02BY

Calibrator blend Expiration Date: 31 JAN 2019

Calibrator blend Lot Number: A0010387

Calibrator blend Catalog Number: C00BZ

Calibrator blend Expiration Date: 30 APR 2018

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

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
Preparer	Xiomara Talavera	Quality Control		06 FEB 2017
Review/Approval	Dawn Stollar	Quality Assurance		06 FEB 2017