



CERTIFICATE OF ANALYSIS

COA - 07377

Product Description: MSD[®] Human Adiponectin Kit

Catalog #/Size: K151BXC-1 / 1 Plate; K151BXC-2 / 5 Plates; K151BXC-3 / 20 Plates

Lot Number: K0013877

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: Z0055484

Plate Lot Catalog Number: N451BXA

Plate Lot Expiration Date: 31 MAR 2018

Antibody blend Lot Number: D0013470

Antibody blend Catalog Number: D21BX

Antibody blend Expiration Date: 31 MAR 2018

Calibrator blend Lot Number: A0010384

Calibrator blend Catalog Number: C01BX

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		05 OCT 2016
Review/Approval	Dawn Stollar	Quality Assurance		06 OCT 2016