



CERTIFICATE OF ANALYSIS

COA - 07381

Product Description: MSD® Human Leptin/Insulin Kit

Catalog #/Size: K15164C-1 / 1 Plate; K15164C-2 / 5 Plates; K15164C-3 / 20 Plates

Lot Number: K0013921

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: Z0055575
Plate Lot Catalog Number: N45164A
Plate Lot Expiration Date: 31 AUG 2018
Antibody blend Lot Number: D0013524
Antibody blend Catalog Number: D21BY
Antibody blend Expiration Date: 31 JUL 2018
Antibody blend Lot Number: D0013525
Antibody blend Catalog Number: D21BZ
Antibody blend Expiration Date: 31 JUL 2018
Calibrator blend Lot Number: A0010393
Calibrator blend Catalog Number: C01BY
Calibrator blend Expiration Date: 30 SEP 2018
Calibrator blend Lot Number: A0010380
Calibrator blend Catalog Number: C00BZ
Calibrator blend Expiration Date: 30 NOV 2017

Specifications:

- Release method: Coating uniformity
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.

Table with 5 columns: Name, Function, Signature, Date. Rows include Preparer (Xiomara Talavera) and Review/Approval (Dawn Stollar).