



CERTIFICATE OF ANALYSIS

COA - 07397

Product Description: MSD® Mouse MCP-1 Ultra-Sensitive Kit
Catalog #/Size: K152AYC-1 / 1 Plate; K152AYC-2 / 5 Plates; K152AYC-4 / 25 Plates
Lot Number: K0035233
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: Z0046192
Plate Lot Catalog Number: N45021B
Plate Lot Expiration Date: 31 JAN 2019
Antibody blend Lot Number: D0033679
Antibody blend Catalog Number: D20AY
Antibody blend Expiration Date: 31 JAN 2019
Antibody source (host/species): Mouse
Calibrator blend Lot Number: A0032256
Calibrator blend Catalog Number: C0021
Calibrator blend Expiration Date: 31 JAN 2019
Calibrator source (host/species): E. coli

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).