



CERTIFICATE OF ANALYSIS

COA - 04439

Product Description: MSD® Human Chemokine 9-Plex Tissue Culture Kit
Catalog #/Size: K15001B-1 / 1 Plate; K15001B-2 / 5 Plates; K15001B-4 / 25 Plates
Lot Number: K0034083
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: Z0044704
Plate Lot Catalog Number: N05001B
Plate Lot Expiration Date: 30 NOV 2015
Antibody blend Lot Number: D0032773
Antibody blend Catalog Number: D2002
Antibody blend Expiration Date: 30 NOV 2015
Antibody source (host/species): Eotaxin, MCP-1, MDC: Mouse
MIP-1β, IP-10, Eotaxin-3, TARC, hIL-8, hMCP-4: Goat
Calibrator blend Lot Number: A0031777
Calibrator blend Catalog Number: C0001
Calibrator blend Expiration Date: 15 Months From Date of Receipt
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 (Rolanda Spratley) and Review/Approval (Dana Short).