



CERTIFICATE OF ANALYSIS

COA - 06284

Product Description: MSD® Human TNF-alpha Ultra-Sensitive Kit
Catalog #/Size: K151BHC-1 / 1 Plate; K151BHC-2 / 5 Plates; K151BHC-4 / 25 Plates
Lot Number: K0035021
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: Z0045882
Plate Lot Catalog Number: L451BHA
Plate Lot Expiration Date: 28 FEB 2018
Antibody blend Lot Number: D0033512
Antibody blend Catalog Number: D21BH
Antibody blend Expiration Date: 28 FEB 2018
Antibody source (host/species): Goat
Calibrator blend Lot Number: A0032176
Calibrator blend Catalog Number: C01BH
Calibrator blend Expiration Date: 18 FEB 2018
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 (Dawn Stollar).