



CERTIFICATE OF ANALYSIS

COA - 06551

Product Description: MSD® Human IL-6R Ultra-Sensitive Kit
Catalog #/Size: K151ALC-1 / 1 Plate; K151ALC-2 / 5 Plates; K151ALC-4 / 25 Plates
Lot Number: K0035081
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: Z0046000
Plate Lot Catalog Number: L451ALA
Plate Lot Expiration Date: 30 JUN 2018
Antibody blend Lot Number: D0033571
Antibody blend Catalog Number: D21AL
Antibody blend Expiration Date: 31 JUL 2018
Antibody source (host/species): Goat
Calibrator blend Lot Number: A0032202
Calibrator blend Catalog Number: C01AL
Calibrator blend Expiration Date: 30 JUN 2018
Calibrator source (host/species): Sf 21

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).