



CERTIFICATE OF ANALYSIS

COA - 07089

Product Description: MSD® Human IL-6 Tissue Culture Kit

Catalog #/Size: K151AKB-1 / 1 Plate; K151AKB-2 / 5 Plates; K151AKB-4 / 25 Plates

Lot Number: K0035160

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: Z0046081
- Plate Lot Catalog Number: L451AKB
- Plate Lot Expiration Date: 30 NOV 2018
- Antibody blend Lot Number: D0033554
- Antibody blend Catalog Number: D21AK
- Antibody blend Expiration Date: 30 APR 2018
- Antibody source (host/species): Goat
- Calibrator blend Lot Number: A0032215
- Calibrator blend Catalog Number: C01XX
- Calibrator blend Expiration Date: 30 JUN 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.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		07 JUL 2016
Review/Approval	Vanessa Castleberry	Quality Assurance		08 JUL 2016