



CERTIFICATE OF ANALYSIS

COA - 06298

Product Description: MSD® Human IL-10 Tissue Culture Kit

Catalog #/Size: K151AOB-1 / 1 Plate; K151AOB-2 / 5 Plates; K151AOB-4 / 25 Plates

Lot Number: K0035052

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: Z0045942
- Plate Lot Catalog Number: N45015B
- Plate Lot Expiration Date: 30 APR 2018
- Antibody blend Lot Number: D0033540
- Antibody blend Catalog Number: D21AO
- Antibody blend Expiration Date: 30 APR 2018
- Antibody source (host/species): Rat
- Calibrator blend Lot Number: A0032192
- Calibrator blend Catalog Number: C0015
- Calibrator blend Expiration Date: 30 APR 2018
- Calibrator source (host/species): GM-CSF: *E. coli*
IL-10, IL-12t/p70: *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.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		10 DEC 2015
Review/Approval	Dawn Stollar	Quality Assurance		15 DEC 2015