



CERTIFICATE OF ANALYSIS

COA - 03968

Product Description: MSD® Human IL-2 Tissue Culture Kit

Catalog #/Size: K151AHB-1 / 1 Plate; K151AHB-2 / 5 Plates; K151AHB-4 / 25 Plates

Lot Number: K0034279

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: Z0044953
- Plate Lot Catalog Number: L451AHB
- Plate Lot Expiration Date: 29 FEB 2016
- Antibody blend Lot Number: D0032936
- Antibody blend Catalog Number: D21AH
- Antibody blend Expiration Date: 29 FEB 2016
- Antibody source (host/species): Mouse
- Calibrator blend Lot Number: A0031874
- Calibrator blend Catalog Number: C01AH
- Calibrator blend Expiration Date: 31 MAR 2016
- 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		26 Nov 2013
Review/Approval	Dana Short	Quality Control		26 Nov 2013