



# CERTIFICATE OF ANALYSIS

COA - 04297

**Product Description:** MSD® Human IFN-gamma Tissue Culture Kit

**Catalog #/Size:** K151AEB-1 / 1 Plate; K151AEB-2 / 5 Plates; K151AEB-4 / 25 Plates

**Lot Number:** K0034475

**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: Z0045173
- Plate Lot Catalog Number: L451AEB
- Plate Lot Expiration Date: 31 JUL 2016
- Antibody blend Lot Number: D0033101
- Antibody blend Catalog Number: D21AE
- Antibody blend Expiration Date: 31 AUG 2016
- Antibody source (host/species): Mouse
- Calibrator blend Lot Number: A0031961
- Calibrator blend Catalog Number: C01AE
- Calibrator blend Expiration Date: 31 AUG 2016
- Calibrator source (host/species): *E. coli*

### Specifications:

- Release method: Coating uniformity
- Mean Intraplate CV:  $\leq 15\%$
- Mean Interplate CV:  $\leq 18\%$
- 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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		03 APR 2014
<b>Review/Approval</b>	Dana Short	Quality Control		18 MAY 2014