



# CERTIFICATE OF ANALYSIS

COA - 06218

**Product Description:** MSD<sup>®</sup> Human IFN-gamma Tissue Culture Kit

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

**Lot Number:** K0034959

**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: Z0045794
- Plate Lot Catalog Number: L451AEB
- Plate Lot Expiration Date: 31 DEC 2017
- Antibody blend Lot Number: D0033464
- Antibody blend Catalog Number: D21AE
- Antibody blend Expiration Date: 31 DEC 2017
- Antibody source (host/species): Mouse
- Calibrator blend Lot Number: A0032149
- Calibrator blend Catalog Number: C01XX
- Calibrator blend Expiration Date: 31 DEC 2017
- 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		18 MAR 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		22 MAR 2016