



# CERTIFICATE OF ANALYSIS

COA - 06989

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

**Catalog #/Size:** K151ADB-1 / 1 Plate; K151ADB-2 / 5 Plates; K151ADB-4 / 25 Plates

**Lot Number:** K0035047

**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: Z0010617

Plate Lot Catalog Number: L15AA

Plate Lot Expiration Date: 31 MAR 2017

Antibody blend Lot Number: D0033535

Antibody blend Catalog Number: D21AD

Antibody blend Expiration Date: 30 APR 2018

Antibody source (host/species): Goat

Antibody blend Lot Number: D0033536

Antibody blend Catalog Number: 3-3009-020001 / 3-3009-020002

Antibody blend Expiration Date: 30 APR 2018

Antibody source (host/species): Goat

Calibrator blend Lot Number: A0032189

Calibrator blend Catalog Number: C01VI

Calibrator blend Expiration Date: 30 APR 2018

Calibrator source (host/species): CHO

### Specifications:

Release method: Coating uniformity

Mean Itraplate 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		07 JUN 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		07 JUN 2016