



## CERTIFICATE OF ANALYSIS

COA - 07642

**Product Description:** MSD® Human TGF-beta1 Kit

**Catalog #/Size:** K151IUC-1 / 1 Plate; K151IUC-2 / 5 Plates; K151IUC-3 / 20 Plates

**Lot Number:** K0050750

**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: Z0055710

Plate Lot Catalog Number: L451IUA

Plate Lot Expiration Date: 28 FEB 2019

Antibody blend Lot Number: D0050687

Antibody blend Catalog Number: D21IU

Antibody blend Expiration Date: 28 FEB 2019

Calibrator blend Lot Number: A0020199

Calibrator blend Catalog Number: N/A

Calibrator blend Expiration Date: 28 FEB 2019

### 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>	Xiomara Talavera	Quality Control		15 DEC 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		16 DEC 2016