



# CERTIFICATE OF ANALYSIS

COA - 06309

**Product Description:** MSD<sup>®</sup> Human TNF-RI Ultra-Sensitive Kit  
**Catalog #/Size:** K151BIC-1 / 1 Plate; K151BIC-2 / 5 Plates; K151BIC-4 / 25 Plates  
**Lot Number:** K0034939  
**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: Z0045751  
 Plate Lot Catalog Number: N45016A  
 Plate Lot Expiration Date: 30 NOV 2017  
 Antibody blend Lot Number: D0033444  
 Antibody blend Catalog Number: D21BI  
 Antibody blend Expiration Date: 30 NOV 2017  
 Antibody source (host/species): Goat  
 Calibrator blend Lot Number: A0032140  
 Calibrator blend Catalog Number: C0016  
 Calibrator blend Expiration Date: 30 NOV 2017  
 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        |
|------------------------|------------------|---------------------|-----------|-------------|
| <b>Preparer</b>        | Rolanda Spratley | Biologics/Chemistry |           | 11 DEC 2015 |
| <b>Review/Approval</b> | Dawn Stollar     | Quality Assurance   |           | 16 DEC 2015 |