



## CERTIFICATE OF ANALYSIS

COA - 07345

**Product Description:** MSD<sup>®</sup> Human RANTES Ultra-Sensitive Kit

**Catalog #/Size:** K151BFC-1 / 1 Plate; K151BFC-2 / 5 Plates; K151BFC-4 / 25 Plates

**Lot Number:** K0035172

**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: Z0046122

Plate Lot Catalog Number: L451BFA

Plate Lot Expiration Date: 31 OCT 2018

Antibody blend Lot Number: D0033639

Antibody blend Catalog Number: D21BF

Antibody blend Expiration Date: 31 OCT 2018

Antibody source (host/species): Goat

Calibrator blend Lot Number: A0032241

Calibrator blend Catalog Number: C01BF

Calibrator blend Expiration Date: 30 SEP 2018

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>	Xiomara Talavera	Quality Control		27 SEP 2016
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance		28 SEP 2016