



## CERTIFICATE OF ANALYSIS

COA - 07060

**Product Description:** MSD<sup>®</sup> Total PRAS40 Whole Cell Lysate Kit

**Catalog #/Size:** K150KRD-1 / 1 Plate; K150KRD-2 / 5 Plates; K150KRD-3 / 20 Plates

**Lot Number:** K0013351

**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: Z0054590

Plate Lot Catalog Number: N450KRA

Plate Lot Expiration Date: 31 OCT 2015

Antibody blend Lot Number: D0012894

Antibody blend Catalog Number: D20KR

Antibody blend Expiration Date: 30 NOV 2015

### 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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		30 JUN 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		30 JUN 2016