



## CERTIFICATE OF ANALYSIS

COA - 06360

**Product Description:** MSD<sup>®</sup> Insulin Signaling Panel (Phospho Protein) Kit

**Catalog #/Size:** K15151C-1 / 1 Plate; K15151C-2 / 5 Plates; K15151C-3 / 20 Plates

**Lot Number:** K0013851

**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: Z0055440

Plate Lot Catalog Number: N45151A

Plate Lot Expiration Date: 31 JAN 2018

Antibody blend Lot Number: D0013439

Antibody blend Catalog Number: D20HP

Antibody blend Expiration Date: 31 JAN 2018

### 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>	Rolanda Spratley	Biologics/Chemistry		03 FEB 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		04 FEB 2016