



## CERTIFICATE OF ANALYSIS

COA - 04408

**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:** K0013568

**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: Z0054943

Plate Lot Catalog Number: N45151A

Plate Lot Expiration Date: 31 AUG 2016

Antibody blend Lot Number: D0013127

Antibody blend Catalog Number: D20HP

Antibody blend Expiration Date: 30 SEP 2016

### 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		08 MAY 2014
<b>Review/Approval</b>	Dana Short	Quality Control		14 MAY 2014