



## CERTIFICATE OF ANALYSIS

COA - 07737

**Product Description:** MSD<sup>®</sup> Phospho (Ser473)/Total Akt Whole Cell Lysate Kit

**Catalog #/Size:** K15100D - Series

**Lot Number:** K0014023

**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: Z0055786

Plate Lot Catalog Number: N45100B

Plate Lot Expiration Date: 31 MAY 2019

Antibody blend Lot Number: D0013520

Antibody blend Catalog Number: D21CB

Antibody blend Expiration Date: 31 JUL 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>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	11 JAN 2017
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 JAN 2017