



# CERTIFICATE OF ANALYSIS

COA - 06727

**Product Description:** MSD<sup>®</sup> Phospho/Total ERK1/2 Whole Cell Lysate Kit  
**Catalog #/Size:** K15107D-1 / 1 Plate; K15107D-2 / 5 Plates; K15107D-3 / 20 Plates  
**Lot Number:** K0013842  
**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: Z0055419  
 Plate Lot Catalog Number: N45107B  
 Plate Lot Expiration Date: 31 JAN 2018  
 Antibody blend Lot Number: D0013430  
 Antibody blend Catalog Number: D21CM  
 Antibody blend Expiration Date: 31 JAN 2018

**Specifications:**

Release method: Coating uniformity  
 Mean Intraplate CV: ≤ 15%  
 Mean Interplate CV: ≤ 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	<i>Rolanda Spratley</i>	01 APR 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	01 APR 2016