



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Akt Signaling (Phospho) Whole Cell Lysate Kit

**Catalog Number:** K15115D - Series

**Lot Number:** K0013952

**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					
<b>Lot #</b>	Z0055638	<b>Catalog #</b>	N45115B	<b>Expiration Date:</b>	30 NOV 2018

Antibody Blend	
<b>Lot Number:</b>	D0013559
<b>Catalog Number:</b>	N/A
<b>Expiration Date:</b>	31 OCT 2018
<b>Storage Temperature:</b>	2-8°C

Antibody Blend	
<b>Lot Number:</b>	D0013560
<b>Catalog Number:</b>	N/A
<b>Expiration Date:</b>	31 OCT 2018
<b>Storage Temperature:</b>	2-8°C

Antibody Blend	
<b>Lot Number:</b>	D0013561
<b>Catalog Number:</b>	N/A
<b>Expiration Date:</b>	31 OCT 2018
<b>Storage Temperature:</b>	2-8°C

Specifications:

Release method:

Coating Uniformity	Specifications
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.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 FEB 2018