



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Phospho(Ser240/244)/Total S6RP Whole Cell Lysate Kit

**Catalog Number:** K15139D - Series

**Lot Number:** K0014144

**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>	Z0056001	<b>Catalog #</b>	N45139A	<b>Expiration Date:</b>	31 DEC 2019

Antibody Blend	
<b>Lot Number:</b>	D0013762
<b>Catalog Number:</b>	D20DH
<b>Expiration Date:</b>	31 DEC 2019
<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>	20 MAR 2018