



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Phospho-4E-BP1(Thr37/46) Whole Cell Lysate Kit

**Catalog Number:** K150KHD - Series

**Lot Number:** K0014225

**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>	Z0056122	<b>Catalog #</b>	N450KHA	<b>Expiration Date:</b>	31 MAY 2020

Antibody Blend	
<b>Lot Number:</b>	D0013850
<b>Catalog Number:</b>	D20KH
<b>Expiration Date:</b>	31 JUL 2020
<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