



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K15115D - Series

Lot Number: K0014290

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 #	Z0056231	Catalog #	N45115B	Expiration Date:	30 SEP 2020

Antibody Blend	
Lot Number:	D0013644
Catalog Number:	N/A
Expiration Date:	30 JUN 2019
Storage Temperature:	2-8°C

Antibody Blend	
Lot Number:	D0013642
Catalog Number:	N/A
Expiration Date:	30 JUN 2019
Storage Temperature:	2-8°C

Antibody Blend	
Lot Number:	D0013923
Catalog Number:	N/A
Expiration Date:	30 SEP 2021
Storage Temperature:	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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	18 SEP 2018