



CERTIFICATE OF ANALYSIS

COA - 07576

Product Description: MSD[®] Phospho/Total ERK1/2 Whole Cell Lysate Kit

Catalog #/Size: K15107D-1 / 1 Plate; K15107D-2 / 5 Plates; K15107D-3 / 20 Plates

Lot Number: K0013993

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: Z0055735

Plate Lot Catalog Number: N45107B

Plate Lot Expiration Date: 31 MAR 2019

Antibody blend Lot Number: D0013601

Antibody blend Catalog Number: N/A

Antibody blend Expiration Date: 31 MAR 2019

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

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
Preparer	Xiomara Talavera	Quality Control		02 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance		12 DEC 2016