



CERTIFICATE OF ANALYSIS

COA - 07298

Product Description: MSD® Phospho(Ser780)/Total Rb Whole Cell Lysate Kit

Catalog #/Size: K15167D-1 / 1 Plate; K15167D-2 / 5 Plates; K15167D-3 / 20 Plates

Lot Number: K0013773

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: Z0055283

Plate Lot Catalog Number: N45167B

Plate Lot Expiration Date: 31 JUL 2017

Antibody blend Lot Number: D0013364

Antibody blend Catalog Number: D20IR

Antibody blend Expiration Date: 31 JUL 2017

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	<i>X Talavera</i>	19 SEP 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	19 SEP 2016