



CERTIFICATE OF ANALYSIS

COA - 05833

Product Description: MSD[®] Phospho(Ser473)/Total Akt Whole Cell Lysate Kit

Catalog #/Size: K15100D-1 / 1 Plate; K15100D-2 / 5 Plates; K15100D-3 / 20 Plates

Lot Number: K0013096

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: Z0054234

Plate Lot Catalog Number: N45100B

Plate Lot Expiration Date: 31 OCT 2013

Antibody blend Lot Number: D0012643

Antibody blend Catalog Number: D21CB

Antibody blend Expiration Date: 31 OCT 2013

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	Rolanda Spratley	Biologics/Chemistry		03 AUG 2015
Review/Approval	Monika Jain	Quality Assurance		03AUG2015