



CERTIFICATE OF ANALYSIS

COA - 03817

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: K0013223

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: Z0054403

Plate Lot Catalog Number: N45100B

Plate Lot Expiration Date: 31 MAR 2014

Antibody blend Lot Number: D0012765

Antibody blend Catalog Number: D21CB

Antibody blend Expiration Date: 30 JUN 2014

Specifications:

Release method: Coating uniformity

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
Preparer	Rolanda Spratley	Biologics/Chemistry		01 AUG 2013
Review/Approval	Dana Short	Quality Control		01 Aug 2013