



CERTIFICATE OF ANALYSIS

COA - 03528

Product Description: MSD[®] Cleaved PARP Whole Cell Lysate Set

Catalog #/Size: C11DE-1

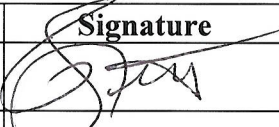

Lot Number: SR627/628

Expiration Date: 28 FEB 2014

Statement: Western blot: This lysate set was tested by western blot using a phospho-specific AKT antibody. Western blot results were consistent with previous lysate preparations. There was no evidence of degraded AKT protein. This lysate set was validated by MSD assay and passed specifications.

Concentration: 2mg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biooigcs/Chemistry		04 MAR 2013
Review/Approval	Michael Eager	Quality Control		20 MAR 2013