



## CERTIFICATE OF ANALYSIS

COA - 04599

**Product Description:** MSD<sup>®</sup> Phospho-Akt (Ser473) Whole Cell Lysate Set

**Catalog #/Size:** C11CA-1



**Lot Number:** CL0010024 / CL0010025

**Expiration Date:** 31 JAN 2018

**Statement:** This lysate set was evaluated by MSD assay and has passed specifications.

Concentration: 2mg/mL

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

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biooigcs/Chemistry		14 AUG 2014
<b>Review/Approval</b>	Dana Short	Quality Control		14 AUG 2014