



## CERTIFICATE OF ANALYSIS

COA – 03547


**Product Description:** Akt Control Pack 1**Catalog Number:** C4115-1**Lot Number:** CP000005**Expiration Date:** 31 OCT 2014**Concentration of Materials Tested:**

	Phospho-Akt (Ser473) (U <sub>P473</sub> /well)	Phospho-Akt (Ser473) (U <sub>P473</sub> /mL)
<b>Akt Control 1</b> (A00C0176)	1.80	72
<b>Akt Control 2</b> (A00C0177)	0.112	4.48

	Total Akt (pg/well)	Total Akt (pg/mL)
<b>Akt Control 1</b> (A00C0176)	485	19 400
<b>Akt Control 2</b> (A00C0177)	21.6	864

<b>Akt Control Pack 1 Components</b>			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Control 1	A00C0176	≤ -70°C	31 OCT 2014
Akt Control 2	A00C0177	≤ -70°C	31 OCT 2014

**Statement:** The above product was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dana Short	Quality Control		14 JAN 2015




# Meso Scale Diagnostics, L.L.C.

Memo: Akt Control Pack 1, Lot CP000005

January 14, 2015

Based on a real-time stability study, the date of expiration for this control pack is extended to 30 months. The expiration date of Lot CP000005 has been extended to April 30, 2015.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		14 Jan 2015

This memo supersedes all other statements of date of expiration for this control pack, Lot CP000005.




# Meso Scale Diagnostics, L.L.C.

Memo: Akt Control 1, Lot A00C0176

January 14, 2015

Based on a real-time stability study, the date of expiration for this control is extended to 30 months. The expiration date of Lot A00C0176 has been extended to April 30, 2015.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		14 Jan 2015

This memo supersedes all other statements of date of expiration for this control, Lot A00C0176.




# Meso Scale Diagnostics, L.L.C.

Memo: Akt Control 2, Lot A00C0177

January 14, 2015

Based on a real-time stability study, the date of expiration for this control is extended to 30 months. The expiration date of Lot A00C0177 has been extended to April 30, 2015.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		14 Jan 2015

This memo supersedes all other statements of date of expiration for this control, Lot A00C0177.