



**RFE MEMO**

Page 1 of 1

**Product Description(s):** Akt Control Pack 1  
**Catalog Number(s) / Size(s):** C4115-1  
**Lot Number:** CP000029

Based on additional testing, the original date of expiration for this control pack has been extended by 48 months. The expiration date of CP000029 has been extended to 31 OCT 2021.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Akt Control 1	A00C0320	31 OCT 2017	31 OCT 2021
Akt Control 2	A00C0321	31 OCT 2017	31 OCT 2021

Function	Name	Signature	Date
Quality	Karen Hamilla		04 NOV 2020

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

COA – 05554

**Product Description:** Akt Control Pack 1**Catalog Number:** C4115-1**Lot Number:** CP000029**Expiration Date:** 31 OCT 2017**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> (A00C0320)	3.77	151
<b>Akt Control 2</b> (A00C0321)	0.204	8.16

	Total Akt (pg/well)	Total Akt (pg/mL)
<b>Akt Control 1</b> (A00C0320)	420	16800
<b>Akt Control 2</b> (A00C0321)	24.6	985

<b>Akt Control Pack 1 Components</b>			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Control 1	A00C0320	≤ -70°C	31 OCT 2017
Akt Control 2	A00C0321	≤ -70°C	31 OCT 2017

**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>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	22 JUL 2016