



**RFE MEMO**

Page 1 of 1

**Product Description(s):** Akt Calibrator  
**Catalog Number(s) / Size(s):** C00MN-2  
**Lot Number:** A0010357

Based on additional testing, the original date of expiration for this calibrator has been extended by 36 months. The expiration date of A0010357 has been extended to 31 Oct 2019.

<b>Function</b>	<b>Name</b>	<b>Signature</b>	<b>Date</b>
<b>Quality Assurance</b>	Dawn Stollar	<i>Dawn Stollar</i>	16 JUL 2018

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



## CERTIFICATE OF ANALYSIS

COA – 06606

**Product Description:** Akt Calibrator

**Catalog Number:** C00MN-2

**Lot Number:** A0010357

**Expiration Date:** 31 OCT 2017

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration of Materials Tested:

	Phospho-Akt (Ser473) (U <sub>P473</sub> /well)	Phospho-Akt (Ser473) (U <sub>P473</sub> /mL)
Calibrator Concentration	872	34 900

	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	87 200	3 490

**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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 MAR 2016