



July 2024

Re: Akt Control Pack 1 (Catalog No.C4115-1), Lot No. CP000110, Revised Concentrations

Dear Valued Customer,

As part of our process for continuous product improvement, we periodically review the performance specifications of MSD products. As we accumulate data from additional lots, if the larger data set indicates a change to a more accurate value, we update the product accordingly.

We recently reviewed the data for the controls in our Akt Control Pack 1, Lot No. CP000110. Based on the larger data set, we decided to reassign the values of the controls for both analytes in this panel. We do not believe that the controls have changed. Rather, we believe that the larger data set provides more accurate values for the controls. The Akt Control Pack 1 is used in the Phospho-Akt (Ser473) Kit and the Total Akt Kit. The revised (Table 1 and Table 2) values and original (Table 3 and Table 4) values for the controls are listed below.

**Table 1:** Phospho-Akt (Ser473) revised concentrations

	<b>Phospho-Akt (Ser473) (UP473/well)</b>	<b>Phospho-Akt (Ser473) (UP473/mL)</b>
<b>AKT Control 1 (A00C0739)</b>	3.27	131
<b>AKT Control 2 (A00C0740)</b>	0.225	9.00

**Table 2:** Total Akt revised concentrations

	<b>Total Akt (pg/well)</b>	<b>Total Akt (pg/mL)</b>
<b>AKT Control 1 (A00C0739)</b>	388	15500
<b>AKT Control 2 (A00C0740)</b>	24.1	964

**Table 3:** Phospho-Akt (Ser473) original concentrations

	<b>Phospho-Akt (Ser473) (UP473/well)</b>	<b>Phospho-Akt (Ser473) (UP473/mL)</b>
<b>AKT Control 1 (A00C0739)</b>	3.65	146
<b>AKT Control 2 (A00C0740)</b>	0.253	10.12



**Table 4:** Total Akt original concentrations

	Total Akt (pg/well)	Total Akt (pg/mL)
<b>AKT Control 1</b> (A00C0739)	405	16200
<b>AKT Control 2</b> (A00C0740)	26	1040

If you have any questions or concerns, please contact MSD Scientific Support at +1 240-314-2798 or [scientificsupport@mesoscale.com](mailto:scientificsupport@mesoscale.com).

Sincerely,

MSD Quality Assurance



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** Akt Control Pack 1


**Catalog Number(s) / Size(s):** C4115-1

**Lot Number:** CP000110

Based on additional testing, the original date of expiration for this control pack has been extended by 24 months. The expiration date of CP000110 has been extended to 30 APR 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Akt Control 1	A00C0739	30 APR 2023	30 APR 2025
Akt Control 2	A00C0740	30 APR 2023	30 APR 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		09 APR 2024

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



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Akt Control Pack

**Catalog Number:** C4115-1

**Lot Number:** CP000110

**Expiration Date:** 30 APR 2023

### Concentration of Materials Tested:

	Phospho-Akt (Ser473) (UP473/well)	Phospho-Akt (Ser473) (UP473/mL)
<b>Akt Control 1</b> (A00C0739)	3.27	130.8
<b>Akt Control 2</b> (A00C0740)	0.225	9

	Total Akt (pg/well)	Total Akt (pg/mL)
<b>Akt Control 1</b> (A00C0739)	387.5	15500
<b>Akt Control 2</b> (A00C0740)	24.1	964

Akt Control Pack 1 Components			
Description	Lot Number	Storage Temperature	Expiration Date
Akt Control 1	A00C0739	≤ -70°C	30 APR 2023
Akt Control 2	A00C0740	≤ -70°C	30 APR 2023

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		10 JUL 2024