



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 1

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)
Akt Control 1 (A00C0739)	3.65	146
Akt Control 2 (A00C0740)	0.253	10.12

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

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
Review/Approval	Karen Hamilla	Quality		08 DEC 2020