



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Akt Control Pack

Catalog Number: C4115-1

Lot Number: CP000193

Expiration Date: 31 DEC 2028

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

Concentration Assignment:

	Phospho-Akt (Ser473) (UP473/well)	Phospho-Akt (Ser473) (UP473/mL)	Total Akt (pg/well)	Total Akt (pg/mL)
Akt Control 1	4.63	185.2	532.5	21300
Akt Control 2	0.275	11.0	32.5	1300

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Akt Control 1	A00C1044	$\leq -70^{\circ}\text{C}$	31 DEC 2028
Akt Control 2	A00C1045	$\leq -70^{\circ}\text{C}$	31 DEC 2028

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 AUG 2026