



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Akt Control Pack

**Catalog Number:** C4115-1

**Lot Number:** CP000174

**Expiration Date:** 31 AUG 2027

### Concentration of Materials Tested:

|                                    | Phospho-Akt (Ser473)<br>(UP473/well) | Phospho-Akt (Ser473)<br>(UP473/mL) |
|------------------------------------|--------------------------------------|------------------------------------|
| <b>Akt Control 1</b><br>(A00C1005) | 2.79                                 | 111.6                              |
| <b>Akt Control 2</b><br>(A00C1006) | 0.196                                | 7.84                               |

|                                    | Total Akt<br>(pg/well) | Total Akt<br>(pg/mL) |
|------------------------------------|------------------------|----------------------|
| <b>Akt Control 1</b><br>(A00C1005) | 465                    | 18600                |
| <b>Akt Control 2</b><br>(A00C1006) | 28.75                  | 1150                 |

| Akt Control Pack 1 Components |            |                            |                 |
|-------------------------------|------------|----------------------------|-----------------|
| Description                   | Lot Number | Storage Temperature        | Expiration Date |
| Akt Control 1                 | A00C1005   | $\leq -70^{\circ}\text{C}$ | 31 AUG 2027     |
| Akt Control 2                 | A00C1006   | $\leq -70^{\circ}\text{C}$ | 31 AUG 2027     |

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  |  | 18 APR 2025 |