



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Akt Calibrator

**Catalog Number:** C00MN-2

**Lot Number:** A0010507


**Expiration Date:** 28 FEB 2025

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

### Concentration Assignment:

	<b>Phospho-Akt (Ser473) (U<sub>P473</sub>/well)</b>	<b>Phospho-Akt (Ser473) (U<sub>P473</sub>/mL)</b>	<b>Total Akt (pg/well)</b>	<b>Total Akt (ng/mL)</b>
Calibrator Concentration	354	14160	110000	4400
Fold Dilution	20X	20X	80X	80X

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

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality		26 SEP 2022