



### CERTIFICATE OF ANALYSIS

**Product Description:** Akt Calibrator  
**Catalog Number:** C00MN-2  
**Lot Number:** A0010521  
**Expiration Date:** 31 JAN 2027  
**Storage Temperature:** ≤ -70°C

	<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	350	14000	107800	4312
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>	Hamilla, Karen	Quality		19 SEP 2024