



CERTIFICATE OF ANALYSIS

Product Description: Akt Calibrator
Catalog Number: C00MN-2
Lot Number: A0010538
Expiration Date: 30 NOV 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

	Phospho-Akt (Ser473) (U _{P473} /well)	Phospho-Akt (Ser473) (U _{P473} /mL)	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	2820	112800	91000	3640
Fold Dilution	20X	20X	80X	80X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 JUL 2026