



CERTIFICATE OF ANALYSIS

Product Description: Akt Calibrator
Catalog Number: C00MN-2
Lot Number: A0010525
Expiration Date: 31 AUG 2027
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	2380	95200	96200	3848
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		18 APR 2025