



CERTIFICATE OF ANALYSIS

Product Description: Akt Calibrator
Catalog Number: C00MN-2
Lot Number: A0010519
Expiration Date: 30 SEP 2026
Storage Temperature: ≤ -70°C

	Phospho-Akt (Ser473) (U_{P473}/well)	Phospho-Akt (Ser473) (U_{P473}/mL)	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	336	13440	110000	4400
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		07 JUN 2024