



CERTIFICATE OF ANALYSIS

Product Description: Akt Calibrator
Catalog Number: C00MN-2
Lot Number: A0010522
Expiration Date: 30 JUN 2027
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	2380	95200	94400	3776
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		28 JAN 2025