



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Akt Calibrator

Catalog Number: C00MN-2

Lot Number: A0010503

Expiration Date: 31 AUG 2024

Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

	Phospho-Akt (Ser473) (U _{P473} /well)	Phospho-Akt (Ser473) (U _{P473} /mL)	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	376	15040	122400	4896
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		31 MAR 2022