



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Akt Calibrator

Catalog Number: C00MN-2

Lot Number: A0010489

Expiration Date: 31 MAY 2023

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	356	14240	114200	4568
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 DEC 2020