



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Akt Calibrator
Catalog Number: C00MN-2
Lot Number: A0010456
Expiration Date: 31 DEC 2020
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration of Materials Tested:

	Phospho-Akt (Ser473) (U _{P473} /well)	Phospho-Akt (Ser473) (U _{P473} /mL)	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	332	13 280	102 000	4 080
Fold Dilution	20X	20X	80X	80X

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 AUG 2018