



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Akt Calibrator

**Catalog Number:** C00MN-2

**Lot Number:** A0010452

**Expiration Date:** 30 NOV 2020

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

### Concentration of Materials Tested:

	Phospho-Akt (Ser473) (U <sub>P473</sub> /well)	Phospho-Akt (Ser473) (U <sub>P473</sub> /mL)	Total Akt (pg/well)	Total Akt (ng/mL)
Calibrator Concentration	316	12 640	102 600	4104
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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	07 AUG 2018