

## ZM-447439

(Catalog No. M75001)

### Alternative Names

N-[4-[[6-Methoxy-7-[3-(4-morpholinyl)propoxy]-4-quinazolinyl]amino]phenyl]benzamide

### Description

ZM-447439 is a selective inhibitor of aurora kinases ARK-1 (A) (IC<sub>50</sub>=1000nM, K<sub>i</sub>=100nM), ARK-2 (B) (IC<sub>50</sub>= 50nM, K<sub>i</sub>= 0.36nM) and ARK-3 (C), that can induce growth arrest, inhibit histone H3 phosphorylation and effectively activate apoptosis. The compound specifically interacts at the ATP-binding site and at the adjacent cleft, found only in aurora kinases. In the presence of Z-M447439, chromosome alignment, segregation, and cytokinesis during mitosis are unsuccessful. Studies show that ZM-447439 can override the spindle assembly checkpoint and accelerate meiosis I in mouse oocytes.

### CAS Number

331771-20-1

### Chemical Formula

C<sub>29</sub>H<sub>31</sub>N<sub>5</sub>O<sub>5</sub>

### Molecular Weight

513.6

### Size

1 mg

### Purity

>98%

### Solubility

Soluble in DMSO (25 mg/mL)

### Storage Temp

-20°C

### Ordering Information:

#### Products

ZM-447439

#### Size

1 mg

#### Cat. No.

M75000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*