

---

## VX-680 (MK-0457, Tozasertib)

(Catalog No. M73000)

---

### Alternative Names

Tozasertib VX 680

### Description

MK 0457 is a potent pan-Aurora kinase inhibitor but favors Aurora A ( $K_i = 0.6$  nM) over Aurora B ( $K_i = 18$  nM) or Aurora C ( $K_i = 4.6$  nM). It shows selectivity against a panel of more than 190 different protein kinases. MK 0457 effectively inhibits proliferation of several different cell lines of clear cell renal carcinoma ( $IC_{50}$ s  $< 10$   $\mu$ M) and blocks the growth of tumors in a rodent model of cancer (80 mg/kg), inhibiting histone H3 phosphorylation and increasing apoptosis. By depleting Aurora activity, MK 0457 disrupts bipolar spindle formation during mitosis, arresting cell cycle progression at the G2/M phase.

### CAS Number

639089-54-6

### Chemical Formula

$C_{23}H_{28}N_8OS$

### Molecular Weight

464.6

### Size

25 mg

### Purity

$\geq 98\%$

### Solubility

DMSO

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
VX-680 (MK-0457, Tozasertib)	25 mg	M73000 -1

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