

## TAME

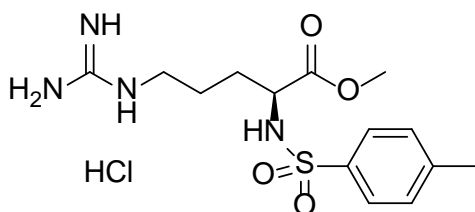
### *N*<sup>2</sup>-[(4-Methylphenyl)sulfonyl]-L-arginine methyl ester hydrochloride

Cat. # SI9840

**Background:** Inhibits the ubiquitin ligase anaphase-promoting complex/cyclosome (APC/C) by inhibiting binding and activation by the APC activators Cdc20 and Cdh1. TAME inhibits cyclin proteolysis in *Xenopus* extracts (IC<sub>50</sub>=12μM) and inhibits Cdc20 association with APC at 200μM. It stabilizes cyclin B1 by prematurely terminating ubiquitination. Substrate for trypsin, thrombin, plasmin, and other proteases.

#### Product Information:

<b>CAS No.</b>	1784-03-8
<b>Purity:</b>	>98% (TLC); NMR (Conforms)
<b>Molecular Weight:</b>	378.9
<b>Physical State:</b>	White Powder
<b>Quantity:</b>	100 mg
<b>Solubility:</b>	May be dissolved in DMSO (35 mg/ml); or water (35 mg/ml)
<b>Storage:</b>	Store desiccated as supplied at ambient temperature for up to 1 year. Store solutions at -20°C for up to 3 months.



**Formula:** C<sub>14</sub>H<sub>22</sub>N<sub>4</sub>O<sub>4</sub>S · HCl

#### References

1. Zeng X. *et al.* Pharmacologic Inhibition of the Anaphase-Promoting Complex Induces A Spindle Checkpoint-Dependent Mitotic Arrest in the Absence of Spindle Damage. *Cancer Cell* 2010 **18**:382
2. Zeng X. *et al.* An APC/C inhibitor stabilizes cyclin B1 by prematurely terminating ubiquitination. *Nat. Chem. Biol.* 2012 **8**:383

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2011 LifeSensors, Inc. All Rights Reserved