

# Triciribine

## Akt inhibitor

Triciribine is a synthetic tricyclic nucleoside which acts as a specific inhibitor of the Akt signaling pathway. It selectively inhibits the phosphorylation and activation of Akt1, -2 and -3 but does not inhibit Akt kinase activity nor known upstream Akt activators such as PI 3-Kinase and PDK1. It inhibits cell growth and induces apoptosis preferentially in cells that express aberrant Akt1. In whole cells triciribine is phosphorylated by adenosine kinase which may be necessary for its activity.

Citations: 7

[View Online »](#)

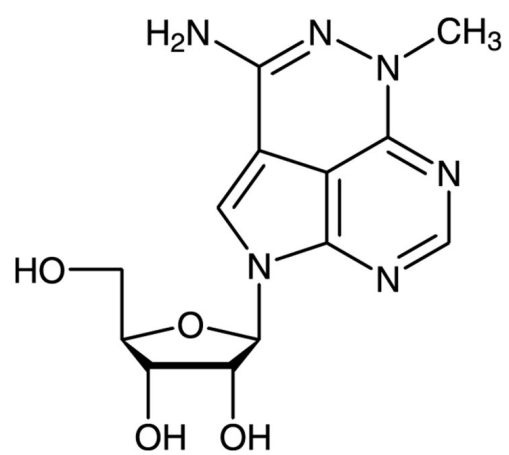
## Ordering Information

[Order Online »](#)

BML-EI332-0001	1mg
BML-EI332-0005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** Pale yellow solid

**CAS** 35943-35-2

**Couple Target** Akt

**Couple Type** Inhibitor

**Formula**  $C_{13}H_{16}N_6O_4$

**MW** 320.3

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO.

**Last modified:** May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)