

Triciribine

Akt inhibitor

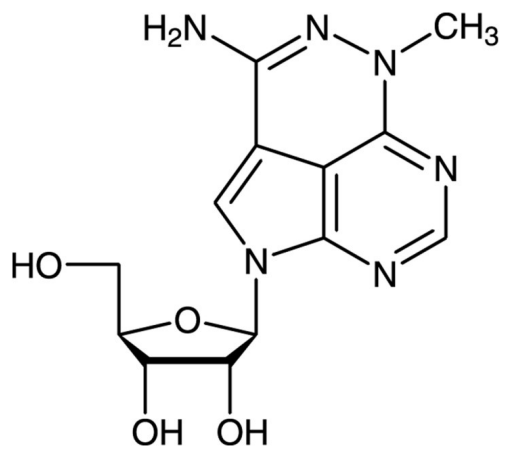
Triciribine is a synthetic tricyclic nucleoside which acts as a specific inhibitor of the Akt signaling pathway. It selectively inhibits the phosphorylation and activation of Akt1, -2 and -3 but does not inhibit Akt kinase activity nor known upstream Akt activators such as PI 3-Kinase and PDK1. It inhibits 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 »](#)

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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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