

# Itraconazole

**CYP 14- $\alpha$  demethylase inhibitor. Autophagy inducer.**

Inhibits 14- $\alpha$  demethylase, a cytochrome P-450 enzyme, thus blocking conversion of lanosterol to ergosterol. Itraconazole inhibits glioblastoma cell proliferation in vitro and in vivo, and stimulates autophagic progression through inhibition of cholesterol trafficking from late endosomes and lysosomes to the plasma membrane. Binds Smoothed (SMO) via a mechanism distinct from that of cyclopamine, inhibiting the Hedgehog pathway. Also inhibits angiogenesis via blocking the binding of VEGF to VEGFR2.

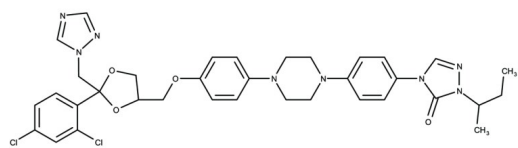
## Ordering Information

[Order Online »](#)

ENZ-CHM265-0250	250mg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored unopened at -20°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Oriconazole, R51211
<b>Appearance</b>	White solid.
<b>CAS</b>	84625-61-6
<b>Couple Target</b>	Cytochrome P450
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{35}H_{36}Cl_2N_8O_4$
<b>Formulation</b>	Lyophilized.
<b>Identity</b>	Identity determined by NMR.
<b>MW</b>	705.6
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in DMSO (1.5 mg/ml, warm).

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-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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