

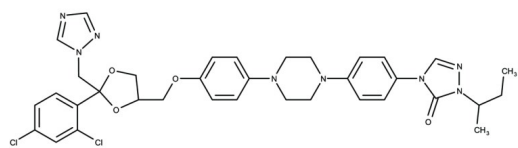
Itraconazole

CYP 14- α demethylase inhibitor. Autophagy inducer.

Inhibits 14- α 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.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

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

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-fr@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