

Indirubin-3'-(2,3-dihydroxypropyl)oximether

Src kinase inhibitor

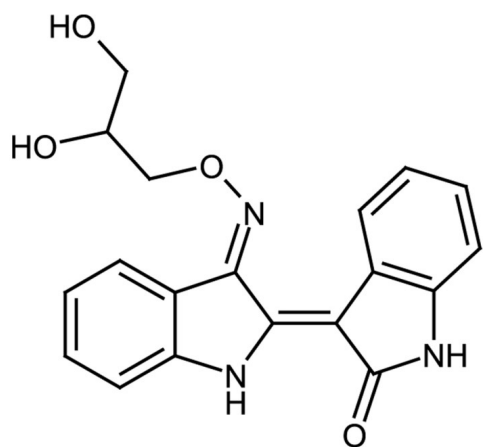
Inhibitor of c-Src ($IC_{50}=0.43\mu\text{M}$) blocking downstream STAT3 signalling. Induces apoptosis in human cancer cells.

Citations: 2

[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

Alternative Name E804

Appearance Dark red solid.

Couple Target Src kinase

Couple Type Inhibitor

Formula $C_{19}H_{17}N_3O_4$

Identity Determined by 1H-NMR and 13C-NMR.

MW 351.4

Purity ≥96% (TLC)

Solubility Soluble in DMSO (1M).

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