

AG-490

JAK inhibitor

AG-490 is a potent inhibitor of the JAK-2 tyrosine kinase. In acute lymphoblastic leukemia (ALL) cells, which abundantly express JAK-2, AG-490 dose-dependently inhibited DNA synthesis, blocked cell growth and induced apoptosis. At 5 μ M, AG-490 almost completely blocked the growth of all pre-B ALL cells but had no significant effect on the growth of mitogen-stimulated normal B or T cells, B-cell lymphoma or T-cell leukemia cells. AG-490 does not significantly inhibit other kinases such as Lck, Lyn, Btk, Syk and Src. It blocks interleukin-7-induced JAK kinase activity in T-cells (JAK-1, JAK-3) and the consequent phosphorylation of PI-3 kinase. AG-490 is cell permeable and is a valuable tool for studying the cellular role of JAK kinases in signal transduction.

Citations: 16

[View Online »](#)

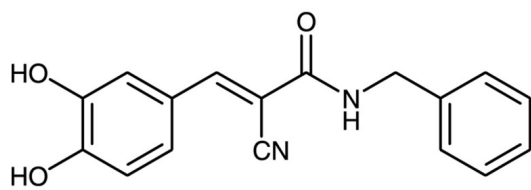
Ordering Information

[Order Online »](#)

BML-EI272-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 2 months.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Benzyl-3,4-dihydroxybenzylidenecyanoacetamide
Appearance	Tan solid.
CAS	133550-30-8
Couple Target	JAK
Couple Type	Inhibitor
Formula	$C_{17}H_{14}N_2O_3$
MW	294.3
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (30mg/ml), dimethyl formamide (40mg/ml) or 100% ethanol (10mg/ml).



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