

SP600125

Jnk inhibitor

Potent, cell permeable, selective, and reversible inhibitor of c-Jun N-terminal kinase (JNK) (IC_{50} =40nM for JNK-1 and JNK-2 and 90nM for JNK-3). The inhibition is competitive with respect to ATP. Exhibits over 300-fold greater selectivity for JNK as compared to ERK1 and p38. Inhibits the phosphorylation of c-Jun and blocks the expression of IL-2, IFN- γ , TNF- α , and COX-2. Blocks IL-1-induced accumulation of phospho-Jun, induction of c-Jun transcription and anti CD3-induced apoptosis of CD4+ CD8+ thymocytes.

Citations: 84

[View Online »](#)

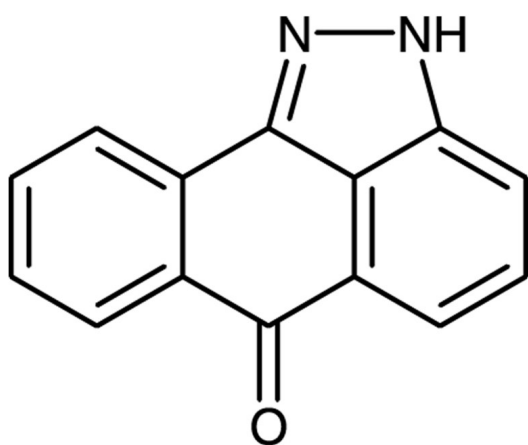
Ordering Information

[Order Online »](#)

BML-EI305-0010	10mg
BML-EI305-0050	50mg

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 at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Anthra(1,9-cd)pyrazol-6(2H)-one, 1,9-Pyrazoloanthrone
Appearance	Yellow to brown solid.
CAS	129-56-6
Couple Target	Jnk
Couple Type	Inhibitor
Formula	$C_{14}H_8N_2O$
MW	220.2
Purity	≥98% (TLC)
RTECS	CB4585000
Solubility	Soluble in DMSO or methanol.
Technical Info / Product Notes	Note: Product is not sterile. Replacement for ADI-HPK-111 .



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