

# 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- $\gamma$ , TNF- $\alpha$ , 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

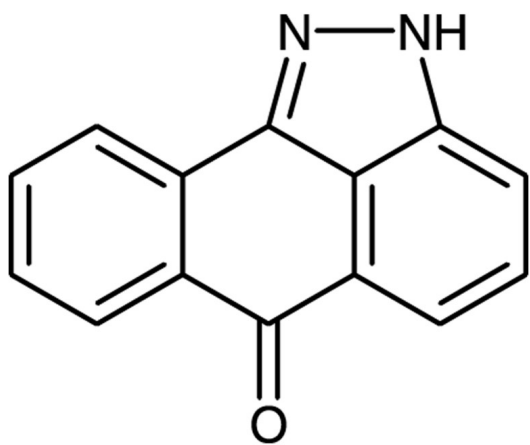
[Order Online »](#)

BML-EI305-0010

10mg

## 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 <b>ADI-HPK-111</b> .



ENZO LIFE SCIENCES,  
Last modified **May 29, 2024**  
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-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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

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

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