

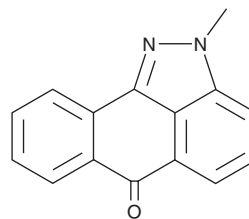
# PRODUCT INFORMATION



## SP600125, negative control

Item No. 18858

**CAS Registry No.:** 54642-23-8  
**Formal Name:** 2-methyl-dibenz[cd,g]indazol-6(2H)-one  
**Synonym:** JNK Inhibitor II, negative control  
**MF:** C<sub>15</sub>H<sub>10</sub>N<sub>2</sub>O  
**FW:** 234.3  
**Purity:** ≥97%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

SP600125, negative control is supplied as a solid. A stock solution may be made by dissolving the SP600125, negative control in the solvent of choice. SP600125, negative control is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

### Description

SP600125 (Item No. 10010466) is a selective, cell-permeable, reversible inhibitor of JNK isoforms 1-3 with K<sub>i</sub> values of 0.19 μM.<sup>1</sup> SP 600125, negative control is a methylated analog of SP600125 that has much lower affinity for JNK isoforms (IC<sub>50</sub>s = 18 and 24 μM for JNK2 and JNK3, respectively).<sup>1</sup>

### Reference

1. Bennett, B.L., Sasaki, D.T., Murray, B.W., *et al.* SP600125, an anthrapyrazolone inhibitor of jun N-terminal kinase. *Proc. Natl. Acad. Sci. USA* **98(24)**, 13681-13686 (2001).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete Safety Data Sheet](#), which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/01/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM