

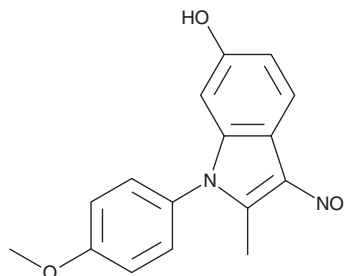
PRODUCT INFORMATION



ID-8

Item No. 15222

CAS Registry No.: 147591-46-6
Formal Name: 1-(4-methoxyphenyl)-2-methyl-3-nitro-1H-indol-6-ol
MF: $C_{16}H_{14}N_2O_4$
FW: 298.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 205, 287, 372 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

ID-8 is supplied as a crystalline solid. A stock solution may be made by dissolving the ID-8 in the solvent of choice, which should be purged with an inert gas. ID-8 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ID-8 in these solvents is approximately 25 and 33.3 mg/ml, respectively.

ID-8 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ID-8 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ID-8 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ID-8 is a cell culture supplement that can sustain self-renewal and pluripotency of mouse embryonic stem cells *in vitro*. At 10 μM it stimulates proliferation in serum-free media at a steady rate for more than 30 days.¹

Reference

1. Miyabayashi, T., Yamamoto, M., Sato, A., *et al.* Indole derivatives sustain embryonic stem cell self-renewal in long-term culture. *Biosci. Biotechnol. Biochem.* **72**(5), 1242-1248 (2008).

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, 11/09/2022

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