

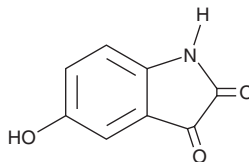
PRODUCT INFORMATION



5-hydroxy Isatin

Item No. 36383

CAS Registry No.: 116569-09-6
Formal Name: 5-hydroxy-1H-indole-2,3-dione
MF: $C_8H_5NO_3$
FW: 163.1
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 214, 255 nm
Supplied as: A 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

5-hydroxy Isatin is supplied as a solid. A stock solution may be made by dissolving the 5-hydroxy isatin in the solvent of choice, which should be purged with an inert gas. 5-hydroxy Isatin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 5-hydroxy isatin in these solvents is approximately 5 and 3 mg/ml, respectively.

Description

5-hydroxy Isatin is an inhibitor of monoamine oxidase A (MAO-A; $IC_{50} = 8.4 \mu\text{M}$).¹ It is selective for MAO-A over MAO-B ($IC_{50} = >100 \mu\text{M}$). 5-hydroxy Isatin ($50 \mu\text{M}$) also inhibits atrial natriuretic peptide-induced activation of particulate guanylyl cyclase in rat brain and heart membranes.²

References

1. Medvedev, A.E., Goodwin, B., Chow, A., *et al.* Inhibitory potency of some isatin analogues on human monoamine oxidase A and B. *Biochem. Pharmacol.* **44**(3), 590-592 (1992).
2. Medvedev, A.E., Goodwin, B.L., Sandler, M., *et al.* Efficacy of isatin analogues as antagonists of rat brain and heart atrial natriuretic peptide receptors coupled to particulate guanylyl cyclase. *Biochem. Pharmacol.* **57**(8), 913-915 (1999).

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, 12/19/2022

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