

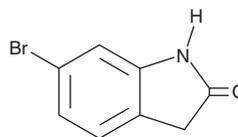
PRODUCT INFORMATION



6-Bromo-2-oxindole

Item No. 20041

CAS Registry No.: 99365-40-9
Formal Name: 6-bromo-1,3-dihydro-2H-indol-2-one
Synonym: 6-Bromooxindole
MF: C₈H₆BrNO
FW: 212.0
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 250 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

6-Bromo-2-oxindole is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-bromo-2-oxindole in the solvent of choice, which should be purged with an inert gas. 6-Bromo-2-oxindole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 6-bromo-2-oxindole in ethanol and DMF is approximately 10 mg/ml and approximately 3 mg/ml in DMSO.

6-Bromo-2-oxindole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 6-bromo-2-oxindole should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 6-Bromo-2-oxindole has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

6-Bromo-2-oxindole is a secondary metabolite that has been isolated from the sea squirt (*D. skoogi*).¹ It is cytotoxic to MDA-MB-231 breast cancer cells with an IC₅₀ value of 74.41 μM. 6-Bromo-2-oxindole has been used in the synthesis of 1,3-disubstituted 2-oxindoles and indolin-2-one p38α inhibitors, which have anti-inflammatory activity.^{2,3}

References

1. Bromley, C.L., Parker-Nance, S., la Mare, J.-A., *et al.* Halogenated oxindole and indoles from the South African marine ascidian *Distaplia skoogi*. *S. Afr. J. Chem.* **66**, 64-68 (2013).
2. Kadin, S.B. 1,3-Disubstituted 2-oxindoles as analgesic and anti-inflammatory agents. **US4,721,712** (1988).
3. Eastwood, P., González, J., González, E., *et al.* Indolin-2-one p38α inhibitors II: Lead optimisation. *Bioorg. Med. Chem. Lett.* **21(18)**, 5270-5273 (2011).

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, 11/11/2022

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