

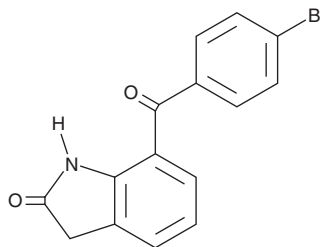
PRODUCT INFORMATION



7-(4-Bromobenzoyl)-1,3-dihydro-2H-indol-2-one

Item No. 43069

CAS Registry No.: 91713-91-6
Synonyms: AHR 10240
MF: $C_{15}H_{10}BrNO_2$
FW: 316.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

7-(4-Bromobenzoyl)-1,3-dihydro-2H-indol-2-one is supplied as a solid. A stock solution may be made by dissolving the 7-(4-bromobenzoyl)-1,3-dihydro-2H-indol-2-one in the solvent of choice, which should be purged with an inert gas. 7-(4-Bromobenzoyl)-1,3-dihydro-2H-indol-2-one is soluble (≥ 10 mg/ml) in DMSO. 7-(4-Bromobenzoyl)-1,3-dihydro-2H-indol-2-one is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

7-(4-Bromobenzoyl)-1,3-dihydro-2H-indol-2-one is a metabolite of bromfenac (Item No. 23763).¹ It is formed from bromfenac via an unstable bromfenac glucuronide intermediate. 7-(4-Bromobenzoyl)-1,3-dihydro-2H-indol-2-one is also a degradation product of bromfenac formed by acid or water hydrolysis, oxidation, and photo or thermal degradation.²

References

1. Driscoll, J.P., Yadav, A.S., and Shah, N.R. Role of glucuronidation and P450 oxidation in the bioactivation of bromfenac. *Chem. Res. Toxicol.* **31**(4), 223-230 (2018).
2. Koppala, S., Reddy, V.R., and Anireddy, J.S. Development and validation of a new stability-indicating RP-UPLC method for the quantitative determination of bromfenac sodium and its impurities in an ophthalmic dosage form. *J. Chromatogr. Sci.* **54**(9) 1514-1521 (2016).

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, 04/07/2025

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