

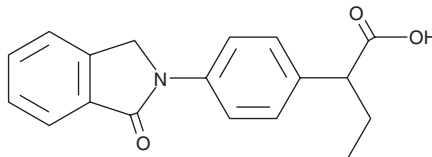
PRODUCT INFORMATION



Indobufen

Item No. 45931

CAS Registry No.: 63610-08-2
Formal Name: 4-(1,3-dihydro-1-oxo-2H-isoindol-2-yl)- α -ethyl-benzeneacetic acid
Synonym: K 3920
MF: $C_{18}H_{17}NO_3$
FW: 295.3
Purity: $\geq 98\%$
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

Indobufen is supplied as a solid. A stock solution may be made by dissolving the indobufen in the solvent of choice, which should be purged with an inert gas. Indobufen is soluble (≥ 10 mg/ml) in DMSO.

Description

Indobufen is a non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor.¹ It reduces thromboxane B_2 (TXB₂) but not prostaglandin E_2 (PGE₂) levels in LPS-stimulated isolated human monocytes. Indobufen inhibits collagen-induced ATP release and platelet aggregation in isolated human platelet-rich plasma (PRP).² It reduces arachidonic acid-induced mortality in mice and rabbits (ED₅₀s = 1.3 and 0.58 mg/kg, respectively).³

References

1. Eligini, S., Violi, F., Banfi, C., *et al.* Indobufen inhibits tissue factor in human monocytes through a thromboxane-mediated mechanism. *Cardiovasc. Res.* **69**(1), 218-226 (2006).
2. Cattaneo, M., Bevilacqua, C., Lecchi, A., *et al.* In vitro and ex vivo effects of indobufen on human platelet aggregation, the release reaction and thromboxane B_2 production. *Haemostasis* **17**(5), 293-300 (1987).
3. Bergamaschi, M., Pierucci, L., and Branzoli, U. Inhibitory activity of indobufen (*) on platelet aggregation in vivo. *Pharmacol. Res. Commun.* **16**(10), 979-985 (1984).

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, 06/19/2026

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