

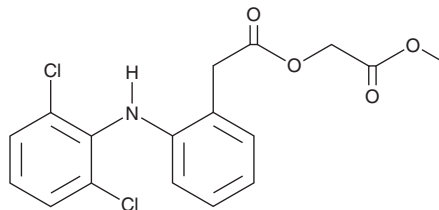
PRODUCT INFORMATION



Aceclofenac methyl ester

Item No. 22121

CAS Registry No.: 139272-66-5
Formal Name: 2-[(2,6-dichlorophenyl)amino]-benzeneacetic acid, 2-methoxy-2-oxoethyl ester
Synonym: Aceclofenac Impurity D
MF: C₁₇H₁₅Cl₂NO₄
FW: 368.2
Purity: ≥98%
UV/Vis.: λ_{max}: 277 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

Aceclofenac methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the aceclofenac methyl ester in the solvent of choice, which should be purged with an inert gas. Aceclofenac methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of aceclofenac methyl ester in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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

Description

Aceclofenac methyl ester is a potential impurity in commercial preparations of the NSAID aceclofenac.¹ It has also been used as a precursor in the synthesis of aceclofenac.²

References

1. Sherikar, O.D., Puranik, M.P., and Yeole, P.G. A validated reversed phase-high performance liquid chromatographic (RP-HPLC) method for simultaneous estimation of aceclofenac drug substance and its related traces impurities in the solid dosage form. *Int. J. Chemtech Res.* **3(2)**, 547-554 (2011).
2. Wu, N.T., Yang, M.S., Wu, H.H., *et al.* New manufacturing method of aceclofenac by deesterification of aceclofenac ester in the presence of metal halide and thiol. *Whail Pharmaceutical Co., Ltd. KR2000009216* (2000).

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, 11/14/2022

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