

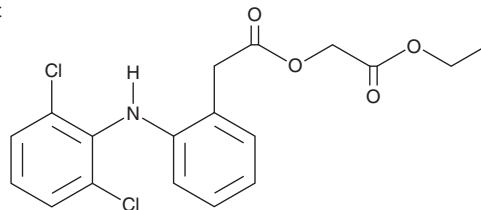
# PRODUCT INFORMATION



## Aceclofenac ethyl ester

Item No. 22115

**CAS Registry No.:** 139272-67-6  
**Formal Name:** 2-[(2,6-dichlorophenyl)amino]-benzeneacetic acid, 2-ethoxy-2-oxoethyl ester  
**MF:** C<sub>18</sub>H<sub>17</sub>Cl<sub>2</sub>NO<sub>4</sub>  
**FW:** 382.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 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 ethyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the aceclofenac ethyl ester in the solvent of choice, which should be purged with an inert gas. Aceclofenac ethyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of aceclofenac ethyl ester in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Aceclofenac ethyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, aceclofenac ethyl ester should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Aceclofenac ethyl 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 ethyl ester is a potential impurity found in commercial preparations of aceclofenac.<sup>1</sup>

### Reference

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).

#### 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