

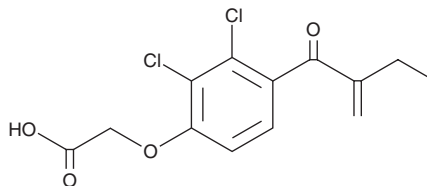
PRODUCT INFORMATION



Ethacrynic Acid

Item No. 26299

CAS Registry No.: 58-54-8
Formal Name: 2-[2,3-dichloro-4-(2-methylene-1-oxobutyl)phenoxy]-acetic acid
Synonyms: MK-595, NSC 85791, NSC 624008
MF: C₁₃H₁₂Cl₂O₄
FW: 303.1
Purity: ≥98%
UV/Vis.: λ_{max}: 268 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Ethacrynic acid (Item No. 26299) is an analytical reference standard categorized as a loop diuretic.¹ Diuretics, including ethacrynic acid, have been used as masking agents in sports doping, including racehorse doping.^{2,3} This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 19536).

References

1. Kunugi, Y., Hiraoka, Y., Hashimoto, Y., *et al.* Ethacrynic acid-sensitive and ATP-dependent Cl⁻ transport in the rat kidney. *Jpn. J. Pharmacol.* **57(2)**, 167-174 (1991).
2. Morra, V., Davit, P., Capra, P., *et al.* Fast gas chromatographic/mass spectrometric determination of diuretics and masking agents in human urine: Development and validation of a productive screening protocol for antidoping analysis. *J. Chromatogr. A* **1135(2)**, 219-229 (2006).
3. Hagedorn, H.W. and Schulz, R. Detection of diuretics in horse urine by GC/MS. *J. Anal. Toxicol.* **16(3)**, 194-198 (1992).

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, 07/17/2023

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