

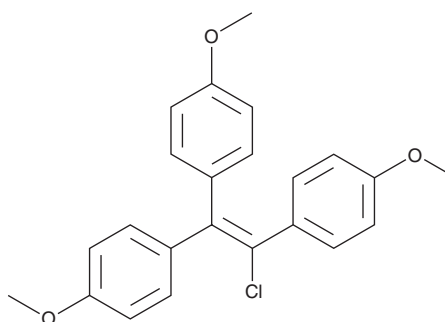
# PRODUCT INFORMATION



## Chlorotrianisene

Item No. 30361

**CAS Registry No.:** 569-57-3  
**Formal Name:** 1,1',1''-(1-chloro-1-ethenyl-2-ylidene)tris[4-methoxy-benzene]  
**Synonyms:** NSC 10108, TACE, Tri-*p*-anisylchloroethylene  
**MF:** C<sub>23</sub>H<sub>21</sub>ClO<sub>3</sub>  
**FW:** 380.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 248, 304 nm  
**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

Chlorotrianisene (TACE) is supplied as a solid. A stock solution may be made by dissolving the TACE in the solvent of choice, which should be purged with an inert gas. TACE is soluble in the organic solvent DMSO.

### Description

TACE is a synthetic non-steroidal estrogen.<sup>1,2</sup> *In vivo*, TACE (5 mg/animal, s.c.) prevents mammary tumor formation induced by demethylbenz[a]anthracene (DMBA) in rats.<sup>1</sup> It also prevents bone loss and uterine atrophy in ovariectomized rats.<sup>2</sup>

### References

1. Kellen, J.A. Suppression of DMBA [7,12-dimethylbenz(a)anthracene] tumor induction in rat mammary glands by pregnancy or chlorotrianisene. *Z. Krebsforsch.* **79(2)**, 75-77 (1973).
2. Zhang, J., Yang, W., and Shen, Z. Prevention of bone loss by chlorotrianisene in oophorectomized rats. *Zhonghua Fu Chan Ke Za Zhi* **32(9)**, 535-537 (1997).

#### 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, 11/16/2022

#### 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