

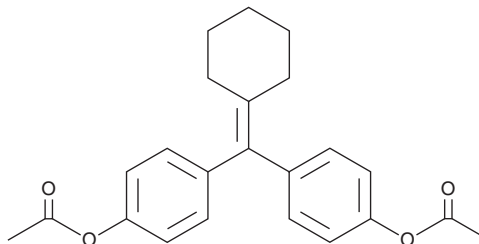
# PRODUCT INFORMATION



## Cyclofenil

Item No. 18676

**CAS Registry No.:** 2624-43-3  
**Formal Name:** 4,4'-(cyclohexylidene)methylene)  
bis-phenol 1,1'-diacetate  
**Synonyms:** ICI 48213, NSC 86464  
**MF:** C<sub>23</sub>H<sub>24</sub>O<sub>4</sub>  
**FW:** 364.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 247 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

Cyclofenil (Item No. 18676) is an analytical reference standard categorized as an anti-estrogen.<sup>1,2</sup> Anti-estrogens, including cyclofenil, have been used as performance-enhancing drugs in sports doping. This product is intended for research and forensic applications.

### References

1. Bowman, S.P., Leake, A. and Morris, I.D. Biological activity and steroid receptor interactions of cyclofenil with the oestrogen target tissues of the brain, pituitary gland and uterus of the rat. *J. Reprod. Fertil.* **65**(2), 355-366 (1982).
2. Gartner, P., Hofbauer, K., Reichel, C., et al. Synthesis and identification of hydroxylated metabolites of the anti-estrogenic agent cyclofenil *J. Mass. Spectrom.* **43**(7), 958-964 (2008).

#### 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, 06/16/2023

#### 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