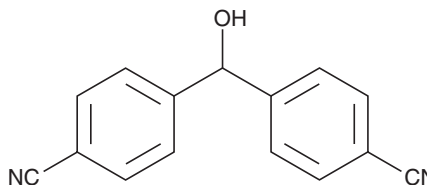


PRODUCT INFORMATION

4,4'-Dicyanobenzhydrol

Item No. 21316

CAS Registry No.: 134521-16-7
Formal Name: 4,4'-(hydroxymethylene)bis-benzonitrile
Synonyms: bis-(4-Cyanophenyl)methanol, bis(p-Cyanophenyl)methanol, CGP 44,645
MF: C₁₅H₁₀N₂O
FW: 234.3
Purity: ≥98%
UV/Vis.: λ_{max}: 242 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

4,4'-Dicyanobenzhydrol (Item No. 21316) is an analytical reference standard that is a metabolite of letrozole (Item No. 11568), an aromatase inhibitor that has been abused during sports doping cycles to reduce gynecomastia associated with androgen abuse.¹ 4,4'-Dicyanobenzhydrol has been used as a marker of letrozole use in athletes. This product is intended for research and forensic applications.

Reference

1. Vari, C.-E., Osz, B.-E., Miklos, A., *et al.* Aromatase inhibitors in men – off-label use, misuse, abuse and doping. *Farmacia* **64**(4), 813-818 (2016).

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, 07/06/2023

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