

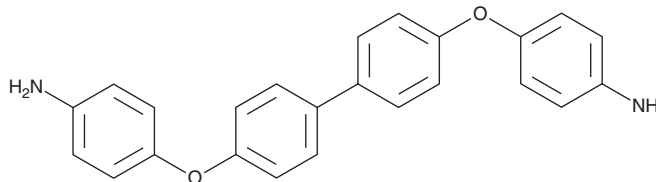
PRODUCT INFORMATION



4,4'-(1,1'-Biphenyl-4,4'-diyldioxy)dianiline

Item No. 39645

CAS Registry No.: 13080-85-8
Formal Name: 4,4'-[[1,1'-biphenyl]-4,4'-diylbis(oxy)]bis-benzenamine
Synonyms: BAPB, 4,4'-Bis(4-aminophenoxy)biphenyl
MF: C₂₄H₂₀N₂O₂
FW: 368.4
Purity: ≥98%
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

4,4'-(1,1'-Biphenyl-4,4'-diyldioxy)dianiline is supplied as a solid. A stock solution may be made by dissolving the 4,4'-(1,1'-biphenyl-4,4'-diyldioxy)dianiline in the solvent of choice, which should be purged with an inert gas. 4,4'-(1,1'-Biphenyl-4,4'-diyldioxy)dianiline is soluble in DMSO. 4,4'-(1,1'-Biphenyl-4,4'-diyldioxy)dianiline is slightly soluble in acetonitrile and THF.

Description

4,4'-(1,1'-Biphenyl-4,4'-diyldioxy)dianiline is a diamine component of polyimide aerogels, super fibers, and thermoplastic shape memory polymers.¹⁻³

References

1. Guo, H., Dewey, O.S., McCorkle, L.S., *et al.* Polyimide aerogels as lightweight dielectric insulators for carbon nanotube cables. *App. Polymer Mater.* **1(7)**, 1680-1688 (2019).
2. Itatani, H., Itatani, T., and Takeichi, T. Polyimide super-fiber from soluble polyimide. *J. Photopolymer Sci. Tech.* **26(3)**, 291-295 (2013).
3. Jayalath, S., Herath, M., Epaarachchi, J., *et al.* Durability and long-term behaviour of shape memory polymers and composites for the space industry - A review of current status and future perspectives. *Polym. Degrad. Stab.* **211**, 110297 (2023).

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, 02/12/2024

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