

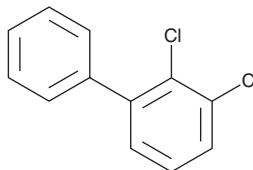
PRODUCT INFORMATION



2,3-Dichlorobiphenyl

Item No. 35886

CAS Registry No.: 16605-91-7
Formal Name: 2,3-dichloro-1,1'-biphenyl
Synonym: PCB 5
MF: $C_{12}H_8Cl_2$
FW: 223.1
Purity: $\geq 98\%$
Supplied as: A solution in methyl acetate
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

2,3-Dichlorobiphenyl is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. 2,3-Dichlorobiphenyl is soluble in chloroform.

Description

2,3-Dichlorobiphenyl is a polychlorobiphenyl (PCB).¹ It decreases the levels of dopamine in PC12 adrenal medulla cells ($EC_{50} = 173 \mu M$). 2,3-Dichlorobiphenyl (10 $\mu g/L$) increases the percentage of *X. laevis* intersex males.² It has been found as a contaminant in urban soil samples.³

References

1. Shain, W., Bush, B., and Seegal, R. Neurotoxicity of polychlorinated biphenyls: Structure-activity relationship of individual congeners. *Toxicol. Appl. Pharmacol.* **111**(1), 33-42 (1991).
2. Qin, Z.-F., Zhou, J.-M., Chu, S.-G., *et al.* Effects of Chinese domestic polychlorinated biphenyls (PCBs) on gonadal differentiation in *Xenopus laevis*. *Environ. Health Perspect.* **111**(4), 553-556 (2003).
3. Cachada, A., Lopes, L.V., Hursthouse, A.S., *et al.* The variability of polychlorinated biphenyls levels in urban soils from five European cities. *Environ. Pollut.* **157**(2), 511-518 (2009).

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, 02/21/2024

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