

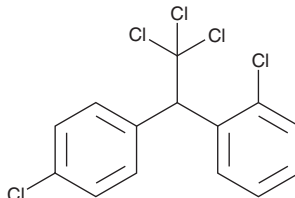
PRODUCT INFORMATION



o,p'-DDT

Item No. 24242

CAS Registry No.: 789-02-6
Formal Name: 1-chloro-2-[2,2,2-trichloro-1-(4-chlorophenyl)ethyl]-benzene
Synonyms: 2,4'-DDT, (\pm)-*o,p'*-DDT, *o,p'*-Dichlorodiphenyltrichloroethane, NSC 33446, NSC 57644
MF: C₁₄H₉Cl₅
FW: 354.5
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 235 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

o,p'-DDT is supplied as a solid. A stock solution may be made by dissolving the *o,p'*-DDT in the solvent of choice. *o,p'*-DDT is slightly soluble in the organic solvent chloroform.

Description

o,p'-DDT is an organochlorine pesticide that induces mortality of malaria mosquito (*A. quadrimaculatus*) fourth-instar larvae and *A. aegypti* larvae when used at concentrations ranging from 0.005 to 0.03 and 0.025 to 5 ppm, respectively.^{1,2} It induces 100% mortality of goldfish (*C. auratus*) at a concentration of 4 ppm.² *o,p'*-DDT is estrogenic, increasing oviduct weight and glycogen content of chicken and Japanese quail.³ It enhances mammary gland differentiation and increases epithelial cell proliferation in mammary terminal end buds in pubertal female rats.⁴ *o,p'*-DDT is a persistent organic pollutant (POP) and is elevated in the sera of pregnant women in malaria-endemic regions of South Africa.⁵

References

1. Jones, H.A., Incho, H.H., and Deonier, C.C. Comparative toxicity of *p,p'*- and *o,p'*-DDT to larvae of *Anopheles quadrimaculatus*. *J. Econ. Entomol.* **39**(5), 672 (1946).
2. Ginsburg, J.M. Comparative toxicity of DDT isomers and related compounds to mosquito larvae and fish. *Science* **105**(2722), 233-234 (1947).
3. Bitman, J., Cecil, H.C., Harris, S.J., *et al.* Estrogenic activity of *o,p'*-DDT in the mammalian uterus and avian oviduct. *Science* **162**(3851), 371-372 (1968).
4. Brown, N.M. and Lamartiniere, C.A. Xenoestrogens alter mammary gland differentiation and cell proliferation in the rat. *Environ. Health Perspect.* **103**(7-8), 708-713 (1995).
5. Röllin, H.B., Sandanger, T.M., Hansen, L., *et al.* Concentration of selected persistent organic pollutants in blood from delivering women in South Africa. *Sci. Total. Environ.* **408**(1), 146-152 (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, 07/12/2023

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