

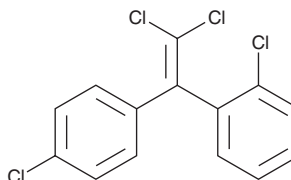
PRODUCT INFORMATION



o,p'-DDE

Item No. 24235

CAS Registry No.: 3424-82-6
Formal Name: 1-chloro-2-[2,2-dichloro-1-(4-chlorophenyl)ethenyl]-benzene
Synonyms: 2,4'-DDE,
o,p'-Dichlorodiphenyldichloroethylene,
NSC 59908
MF: C₁₄H₈Cl₄
FW: 318.0
Purity: ≥95%
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'-DDE is supplied as a solid. A stock solution may be made by dissolving the *o,p'*-DDE in the solvent of choice, which should be purged with an inert gas. *o,p'*-DDE is slightly soluble in warm chloroform and methanol.

Description

o,p'-DDE is a metabolite and degradation product of the organochlorine pesticide DDT.¹ It accumulates in smallmouth buffalo, channel catfish, and largemouth bass as well as sediment in the Huntsville Spring Branch-Indian Creek tributary system surrounding a DDT manufacturing plant where it is considered a persistent organic pollutant (POP). *o,p'*-DDE inhibits estrogen binding to rainbow trout estrogen receptors (rtERs) with an IC₅₀ value of 3.2 μM.² It induces concentration-dependent secretion of estradiol by granulosa and theca cell co-cultures isolated from porcine ovarian follicles.³ *In ovo* exposure to *o,p'*-DDE increases degeneration of ovarian follicles and reduces testicular size in Japanese medaka (*O. latipes*).⁴

References

1. Garrison, A.W., Cyterski, M., Roberts, K.D., *et al.* Occurrences and fate of DDT principal isomers/metabolites, DDA, and *o,p'*-DDD enantiomers in fish, sediment and water at a DDT-impacted Superfund site. *Environ. Pollut.* **194**, 224-234 (2014).
2. Matthews, J., Celius, T., Halgren, R., *et al.* Differential estrogen receptor binding of estrogenic substances: A species comparison. *J. Steroid. Biochem. Mol. Biol.* **74(4)**, 223-234 (2000).
3. Wójtowicz, A.K., Gregoraszczyk, E.L., Ptak, A., *et al.* Effect of single and repeated *in vitro* exposure of ovarian follicles to *o,p'*-DDT and *p,p'*-DDT and their metabolites. *Pol. J. Pharmacol.* **56(4)**, 465-472 (2004).
4. Papoulias, D.M., Villalobos, S.A., Meadows, J., *et al.* *In ovo* exposure to *o,p'*-DDE affects sexual development but not sexual differentiation in Japanese medaka (*Oryzias latipes*). *Environ. Health Perspect.* **111(1)**, 29-32 (2003).

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, 11/07/2022

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