

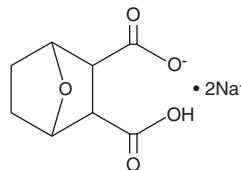
PRODUCT INFORMATION



Endothall (sodium salt)

Item No. 44847

CAS Registry No.: 129-67-9
Formal Name: 7-oxabicyclo[2.2.1]heptane-2,3-dicarboxylic acid, disodium salt
MF: $C_8H_9O_5 \cdot 2Na$
FW: 230.1
Purity: $\geq 95\%$
Supplied as: A solid
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

Endothall (sodium salt) is supplied as a solid. Aqueous solutions of endothall (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. Endothall (sodium salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Endothall is an aquatic plant growth regulator.¹ It reduces curly leaf pondweed appearance frequency in and dry biomass from lakes when applied annually at a concentration of 1 mg/L. Endothall reduces leaf chlorophyll content in *A. thaliana* plants and root length in *A. thaliana* seedlings (IC_{50} s = 1.36 mM and 13.9 μ M, respectively) and decreases duckweed growth rate (IC_{50} = 10 μ M).² It induces production of ethylene in bean leaves when applied at a concentration of 1 μ M.³ Endothall is also an inhibitor of protein phosphatase 2A (PP2A; IC_{50} = 19-50 nM).⁴ It is toxic to mice (LD_{50} = 5-50 mg/kg).

References

1. Johnson, J.A., Jones, A.R., and Newman, R.M. Evaluation of lakewide, early season herbicide treatments for controlling invasive curlyleaf pondweed (*Potamogeton crispus*) in Minnesota lakes. *Lake Reserv. Manag.* **28**(4), 346-363 (2012).
2. Bajsa, J., Pan, Z., Dayan, F.E., et al. Validation of serine/threonine protein phosphatase as the herbicide target site of endothall. *Pestic. Biochem. Physiol.* **102**(1), 38-44 (2012).
3. Abeles, A.L. and Abeles, F.B. Biochemical pathway of stress-induced ethylene. *Plant Physiol.* **50**(4), 496-498 (1972).
4. Li, Y.-M. and Casida, J.E. Cantharidin-binding protein: Identification as protein phosphatase 2A. *Proc. Natl. Acad. Sci. USA* **89**(24), 11867-11870 (1992).

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, 04/14/2026

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