

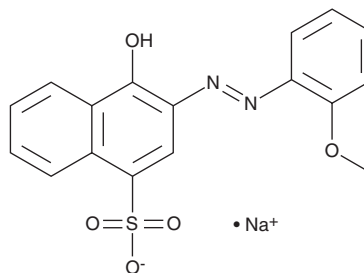
PRODUCT INFORMATION



Acid Red 4

Item No. 44621

CAS Registry No.: 5858-39-9
Formal Name: 4-hydroxy-3-[2-(2-methoxyphenyl) diazenyl]-1-naphthalenesulfonic acid, monosodium salt
Synonyms: Azoeosin C.I. No. 14710, Amacid Eosine, Azo Eosin, Azo Eosine G, C.I. 14710
MF: $C_{17}H_{13}N_2O_5S \cdot Na$
FW: 380.4
UV/Vis.: λ_{max} : 240, 306, 376, 506 nm
Abs. Max.: 508 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

Acid red 4 is supplied as a solid. A stock solution may be made by dissolving the acid red 4 in the solvent of choice, which should be purged with an inert gas. Acid red 4 is soluble (≥ 10 mg/ml) in DMSO.

Description

Acid red 4 is an azo dye.¹ It displays an absorbance maximum at 508 nm. Formulations containing acid red 4 have been used as textile dyes.

Reference

1. Coughlin, M.F., Kinkle, B.K., and Bishop, P.L. Degradation of azo dyes containing aminonaphthol by *Sphingomonas* sp strain 1CX. *J. Ind. Microbiol. Biotechnol.* **23**(4-5), 341-346 (1999).

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, 12/03/2025

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