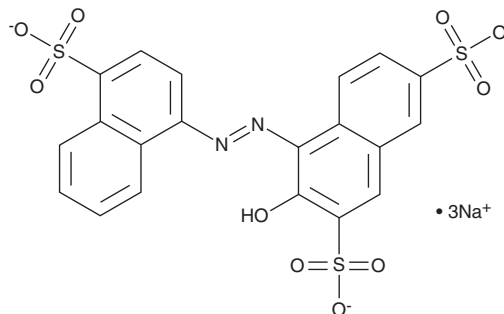


PRODUCT INFORMATION



Acid Red 27 Item No. 44512

CAS Registry No.: 915-67-3
Formal Name: 3-hydroxy-4-[2-(4-sulfo-1-naphthalenyl) diazenyl]-2,7-naphthalenedisulfonic acid, trisodium salt
Synonyms: Amaranth, AR27, C.I. 16185, C.I. Acid Red 27, FD&C Red No. 2
MF: $C_{20}H_{11}N_2O_{10}S_3 \cdot 3Na$
FW: 538.5
UV/Vis.: λ_{max} : 218, 524 nm
Abs. Max.: 520 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 27 is supplied as a solid. Aqueous solutions of acid red 27 can be prepared by directly dissolving the solid in aqueous buffers. Acid red 27 is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Acid red 27 is an azo dye.¹ It displays an absorption maximum at 520 nm. Acid red 27 (47 mg/kg) increases total lymphocytes and basophils and decreases total neutrophils and monocytes but does not affect body weight or serum protein levels in rats.² It has been found as a contaminant in wastewater.¹ Formulations containing acid red 27 have previously been used as food and cosmetic dyes.

References

1. Lehl, H.K., Ong, S.-A., Ho, L.-N., *et al.* Decolourization and mineralization of Acid Red 27 metabolites by using multiple zoned aerobic and anaerobic constructed wetland reactor. *Desalination Water Treat.* **160**, 81-93 (2019).
2. Hashem, M.M., Atta, A.H., Arbid, M.S., *et al.* Immunological studies on Amaranth, Sunset Yellow and Curcumin as food colouring agents in albino rats. *Food Chem. Toxicol.* **48(6)**, 1581-1586 (2010).

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, 12/16/2025

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