

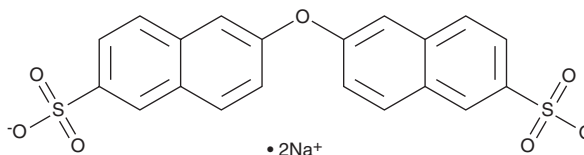
PRODUCT INFORMATION



6,6'-Oxybis(2-naphthalenesulfonate) (sodium salt)

Item No. 45595

CAS Registry No.: 61551-82-4
Formal Name: 6,6'-oxybis-2-naphthalenesulfonic acid, disodium salt
Synonym: DONS
MF: C₂₀H₁₂O₇S₂ • 2Na
FW: 474.4
Purity: ≥98%
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.

Description

6,6'-Oxybis(2-naphthalenesulfonate) is a byproduct in the synthesis of FD&C Yellow No. 6 and FD&C Red No. 40 (Allura Red AC; Item No. 44514).^{1,2} It is formed during the synthesis of Schaeffer's salt, an intermediate used in the synthesis of FD&C Yellow No. 6.¹ 6,6'-Oxybis(2-naphthalenesulfonate) is also an impurity formed in the synthesis of FD&C Yellow No. 5.³

References

1. Bailey, J.E. and Calvey, R.J. Spectral compilation of dyes, intermediates, and other reaction products structurally related to FD&C Yellow No. 6. *J. Assoc. Off. Anal. Chem.* **58(6)**, 1087-1128 (2020).
2. Cox, E.A. and Reed, G.F. High performance liquid chromatographic determination of intermediates and two reaction by-products in FD&C Red No. 40: Collaborative study. *J. Assoc. Off. Anal. Chem.* **64(2)**, 324-331 (2020).
3. Ogura, N., Nakamura, Y., Shinomiya, K., et al. Separation of impurities in commercial food Yellow No. 5 by counter alternative current chromatography and structural analyses. *Anal. Sci.* **11**, 759-763 (1994).

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, 05/12/2026

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