

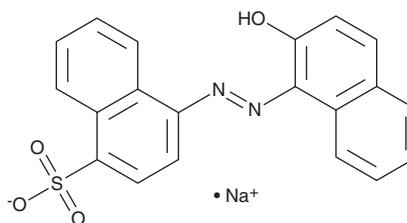
PRODUCT INFORMATION



Acid Red 88

Item No. 44524

CAS Registry No.: 1658-56-6
Formal Name: 4-[2-(2-hydroxy-1-naphthalenyl) diazenyl]-1-naphthalenesulfonic acid, monosodium salt
Synonyms: AR88, C.I. 15620, Rocceline
MF: $C_{20}H_{13}N_2O_4S \cdot Na$
FW: 400.4
Absorbance: 505 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 88 is supplied as a solid. A stock solution may be made by dissolving the acid red 88 in the solvent of choice, which should be purged with an inert gas. Acid red 88 is soluble (≥10 mg/ml) in DMSO.

Description

Acid red 88 is an azo dye.¹ It displays an absorbance maximum at 505 nm. Acid red 88 has been found as a contaminant in textile wastewaters. Formulations containing acid red 88 have been used as textile dyes.

Reference

1. Song, Y.-L., Li, J.-T., and Chen, H. Degradation of C.I. Acid Red 88 aqueous solution by combination of Fenton's reagent and ultrasound irradiation. *J. Chem. Technol. Biotechnol.* **84**(4), 578-583 (2008).

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, 12/17/2025

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