

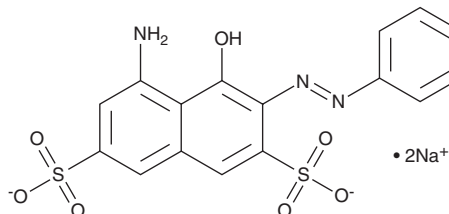
PRODUCT INFORMATION



Acid Red 33

Item No. 44709

CAS Registry No.: 3567-66-6
Formal Name: 5-amino-4-hydroxy-3-(2-phenyldiazenyl)-2,7-naphthalenedisulfonic acid, disodium salt
Synonyms: AR33, D&C Red No. 33
MF: $C_{16}H_{11}N_3O_7S_2 \cdot 2Na$
FW: 467.4
Abs Max: 530 nm
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

Acid red 33 is supplied as a solid. A stock solution may be made by dissolving the acid red 33 in the solvent of choice, which should be purged with an inert gas. Acid red 33 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of acid red 33 can be prepared by directly dissolving the solid in aqueous buffers. Acid red 33 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

Acid red 33 is an azo dye.¹ It displays an absorbance maximum at 530 nm. Formulations containing acid red 33 have been used in dental care products and cosmetics.

Reference

1. Nabizadeh Chianeh, F., Basiri Parsa, J., and Rezaei Vahidian, H. Artificial neural network modeling for removal of azo dye from aqueous solutions by Ti anode coated with multiwall carbon nanotubes. *Environ. Prog. Sustainable Energy* **36**(6), 1778-1784 (2017).

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, 11/06/2025

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