

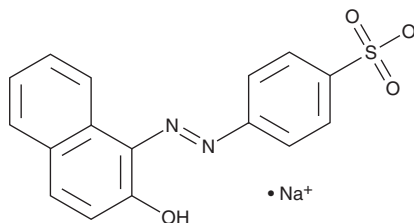
PRODUCT INFORMATION



Acid Orange 7

Item No. 44523

CAS Registry No.: 633-96-5
Formal Name: 4-[2-(2-hydroxy-1-naphthalenyl)diazenyl]-benzenesulfonic acid, monosodium salt
Synonyms: AO7, Orange 4, Orange II, 2-Naphthol Orange
MF: $C_{16}H_{11}N_2O_4S \cdot Na$
FW: 350.3
Abs./Em. Max.: 480/580 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 orange 7 is supplied as a solid. A stock solution may be made by dissolving the acid orange 7 in the solvent of choice, which should be purged with an inert gas. Acid orange 7 is soluble (≥10 mg/ml) in DMSO.

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 orange 7 can be prepared by directly dissolving the solid in aqueous buffers. Acid orange 7 is sparingly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Acid orange 7 is an azo dye.¹ It displays absorbance/emission maxima of 480/580 nm, respectively. Formulations containing acid orange 7 have been used as dyes in fabric, food, and cosmetics.

Reference

1. Vinodgopal, K. and Kamat, P.V. Photochemistry of textile azo dyes. Spectral characterization of excited state, reduced and oxidized forms of Acid Orange 7. *J. Photochem. Photobiol. A Chem.* **83**(2), (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, 10/30/2025

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