

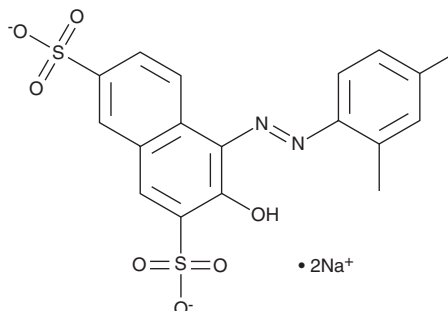
PRODUCT INFORMATION



Acid Red 26

Item No. 44531

CAS Registry No.: 3761-53-3
Formal Name: 4-[2-(2,4-dimethylphenyl)diazenyl]-3-hydroxy-2,7-naphthalenedisulfonic acid, disodium salt
Synonyms: Acid Scarlet GR, AR-26, Ponceau MX, Ponceau Xylidine, Xylidine Ponceau 2R
MF: $C_{18}H_{14}N_2O_7S_2 \cdot 2Na$
FW: 480.4
Abs. Max: 500 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 26 is supplied as a solid. Aqueous solutions of acid red 26 can be prepared by directly dissolving the solid in aqueous buffers. Acid red 26 is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Acid red 26 is an azo dye.¹ It displays an absorbance maximum at 500 nm. Acid red 26 induces mortality in zebrafish larvae at 2,800 µg/ml and decreases heart rate and induces cardiac edema, yolk sac absorption delay, and swimming bladder deflation in zebrafish larvae in a concentration-dependent manner.² It increases sister chromatid exchange (SCE) in mouse bone marrow cells when administered at doses of 62.5 and 125 mg/kg.³ Formulations containing acid red 26 have been used as textile dyes.

References

1. Ma, Q., Bai, H., Zhang, Q., *et al.* Determination of carcinogenic and allergenic dyestuffs in toys by LC coupled to UV/Vis spectrometry and tandem mass spectrometry. *Chromatographia* **72**(1), 85-93 (2010).
2. Shen, B., Liu, H.-C., Ou, W.-B., *et al.* Toxicity induced by basic violet 14, direct red 28 and acid red 26 in zebrafish larvae. *J. Appl. Toxicol.* **35**(12), 1473-1480 (2015).
3. Parodi, S., Zunino, A., Ottaggio, L., *et al.* Lack of correlation between the capability of inducing sister-chromatid exchanges in vivo and carcinogenic potency, for 16 aromatic amines and azo derivatives. *Mutat. Res.* **108**(1-3), 225-238 (1983).

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, 11/10/2025

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