

PRODUCT INFORMATION



Direct Red 80

Item No. 39535

CAS Registry No.: 2610-10-8
Formal Name: 7,7'-(carbonyldiimino)bis[4-hydroxy-3-[2-[2-sulfo-4-[2-(4-sulphophenyl)phenyl]diazenyl]-2-naphthalenesulfonic acid, hexasodium salt

Synonyms: C.I. 35780, Sirius Red, Sirius Red F3B

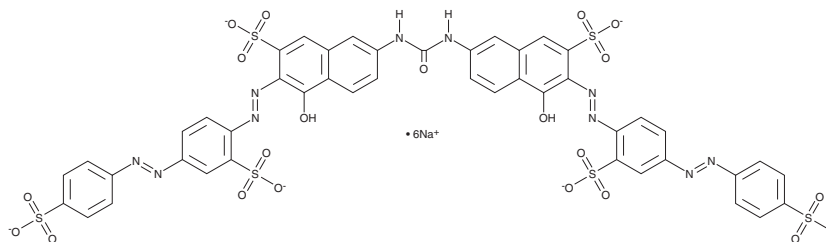
MF: $C_{45}H_{26}N_{10}O_{21}S_6 \cdot 6Na$
FW: 1,373.1

UV/Vis.: λ_{max} : 280, 528 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

Direct red 80 is supplied as a solid. A stock solution may be made by dissolving the direct red 80 in the solvent of choice, which should be purged with an inert gas. Direct red 80 is slightly soluble in DMSO and dimethyl formamide.

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 direct red 80 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of direct red 80 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Direct red 80 is an anionic hexasulfonic acid dye.¹ It displays an absorbance maximum of 540 nm. Direct red 80 has been used as a crosslinking dye for collagen or amyloids, as well as a dye for fabrics.^{1,2}

References

1. Junqueira, L.C., Bignolas, G., and Brentani, R.R. Picrosirius staining plus polarization microscopy, a specific method for collagen detection in tissue sections. *Histochem. J.* **11**(4), 447-455 (1979).
2. Dapson, R.W., Fagan, C., Kiernan, J.A., et al. Certification procedures for sirius red F3B (CI 35780, direct red 80). *Biotech. Histochem.* **86**(3), 133-139 (2011).

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, 07/02/2024

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