

PRODUCT INFORMATION

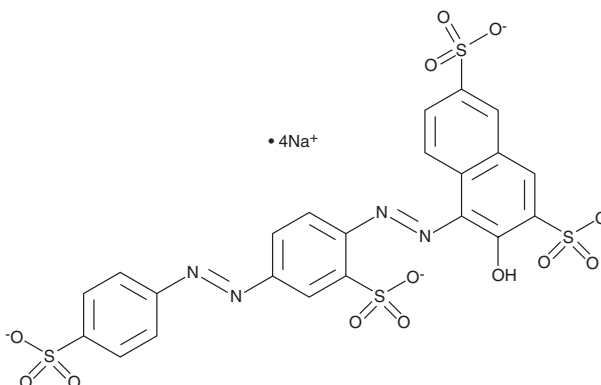


Ponceau S (sodium salt)

Item No. 14330

CAS Registry No.: 6226-79-5
Formal Name: 3-hydroxy-4-[2-[2-sulfo-4-[2-(4-sulfophenyl)diazenyl]phenyl]diazenyl]-2,7-naphthalenedisulfonic acid, tetrasodium salt

Synonym: Acid Red 112
MF: C₂₂H₁₂N₄O₁₃S₄ • 4Na
FW: 760.6
UV/Vis.: λ_{max}: 206, 239, 315, 357, 530 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ponceau S (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the ponceau S (sodium salt) in the solvent of choice, which should be purged with an inert gas. Ponceau S (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ponceau S (sodium salt) in ethanol and DMF is approximately 0.5 mg/ml and approximately 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 ponceau S (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ponceau S (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ponceau S is a stain that permits the rapid, reversible detection of proteins immobilized on various matrices, including nitrocellulose and polyvinylidene fluoride membranes.¹ It can be used to locate proteins or evaluate efficiency of transfer. Reversible staining with ponceau S has also been validated as an alternative to actin probing as a loading control.² This stain can also be used as a fiducial marker in mass spectrometric imaging.³

References

1. Yonan, C.R., Duong, P.T., and Chang, F.N. High-efficiency staining of proteins on different blot membranes. *Anal. Biochem.* **338**(1), 159-161 (2005).
2. Romero-Calvo, I., Ocón, B., Martínez-Moya, P., *et al.* Reversible Ponceau staining as a loading control alternative to actin in Western blots. *Anal. Biochem.* **401**(2), 318-320 (2010).
3. Chughtai, K., Jiang, L., Greenwood, T.R., *et al.* Fiducial markers for combined 3-dimensional mass spectrometric and optical tissue imaging. *Anal. Chem.* **84**(4), 1817-1823 (2012).

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, 12/16/2022

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