

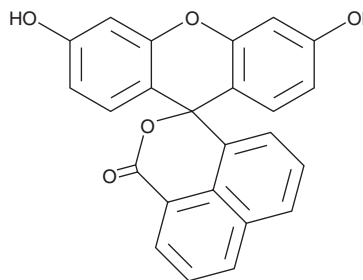
PRODUCT INFORMATION



Resorcinolnaphthalein

Item No. 13082

CAS Registry No.: 41307-63-5
Formal Name: 3',6'-dihydroxy-spiro[1H,3H-naphtho[1,8-cd]pyran-1,9'-[9H]xanthen-3-one
Synonym: NSC 354317
MF: C₂₄H₁₄O₅
FW: 382.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 278 nm
Supplied as: A crystalline 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

Resorcinolnaphthalein is supplied as a crystalline solid. A stock solution may be made by dissolving the resorcinolnaphthalein in the solvent of choice, which should be purged with an inert gas. Resorcinolnaphthalein is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of resorcinolnaphthalein in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Resorcinolnaphthalein is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, resorcinolnaphthalein should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Resorcinolnaphthalein has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Angiotensin-converting enzyme 2 (ACE2) is an enzyme that is cardioprotective and renoprotective. Resorcinolnaphthalein increases the activity of ACE2 *in vitro* (EC₅₀ = 19.5 μM).¹ Activation of ACE2 in rats produces a decrease in blood pressure and improvement in cardiac function. In spontaneously hypertensive rats it produces a reversal of myocardial, perivascular, and renal fibrosis.¹

Reference

1. Prada, J.A.H., Ferreira, A.J., Katovich, M.J., *et al.* Structure-based identification of small-molecule angiotensin-converting enzyme 2 activators as novel antihypertensive agents. *Hypertension* **51**, 1312-1317 (2008).

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, 12/06/2022

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