

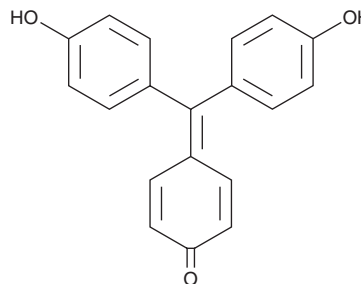
PRODUCT INFORMATION



p-Rosolic Acid

Item No. 44935

CAS Registry No.: 603-45-2
Formal Name: 4-[bis(4-hydroxyphenyl)methylene]-2,5-cyclohexadien-1-one
Synonyms: Aurin, *para*-Rosolic Acid
MF: C₁₉H₁₄O₃
FW: 290.3
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



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

Laboratory Procedures

p-Rosolic acid is supplied as a solid. A stock solution may be made by dissolving the *p*-rosolic acid in the solvent of choice, which should be purged with an inert gas. *p*-Rosolic acid is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

p-Rosolic acid is an acid-base indicator.¹ It is yellow in color below and red in color above its transition range between pH 6.6 and 8.0.

Reference

1. Shokrollahi, A., Gohari, M., and Ebrahimi, F. Determination of acidity constants of *p*-rosolic acid and bromoxynol blue by solution scanometric method. *Anal. Bioanal. Chem. Res.* **5**(1), 67-79 (2018).

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, 02/06/2026

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