

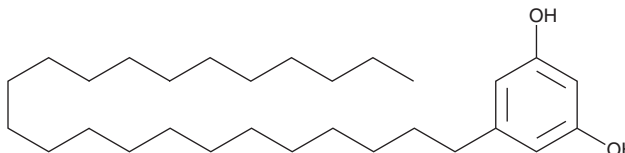
PRODUCT INFORMATION



5-Tricosylresorcinol

Item No. 30032

CAS Registry No.: 70110-60-0
Formal Name: 5-tricosyl-1,3-benzenediol
MF: $C_{29}H_{52}O_2$
FW: 432.7
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years
Item Origin: Plant/Bran powder



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

Laboratory Procedures

5-Tricosylresorcinol is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-tricosylresorcinol in the solvent of choice, which should be purged with an inert gas. 5-Tricosylresorcinol is soluble in ethanol.

Description

5-Tricosylresorcinol is an alkylresorcinol that has been found in wheat, rye, triticale, and barley.¹ It reduces hydrogen peroxide-induced DNA damage in HT-29 cells when used at a concentration of $100 \mu M$.²

References

1. Ross, A.B., Shepherd, M.J., Schüpphaus, M., *et al.* Alkylresorcinols in cereals and cereal products. *J. Agric. Food Chem.* **51**(14), 4111-4118 (2003).
2. Parikka, K., Rowland, I.R., Welch, R.W., *et al.* In vitro antioxidant activity and antigenotoxicity of 5-n-alkylresorcinols. *J. Agric. Food Chem.* **54**(5), 1646-1650 (2006).

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, 01/31/2020

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