

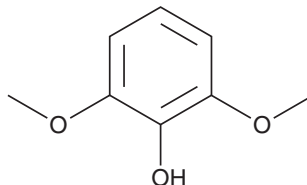
# PRODUCT INFORMATION



## Syringol

Item No. 34475

CAS Registry No.: 91-10-1  
Formal Name: 2,6-dimethoxy-phenol  
Synonyms: 2,6-Dimethoxyphenol,  
Pyrogallol 1,3-dimethyl ether  
MF:  $C_8H_{10}O_3$   
FW: 154.2  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 204, 206 nm  
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Syringol is supplied as a solid. A stock solution may be made by dissolving the syringol in the solvent of choice, which should be purged with an inert gas. Syringol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of syringol in these solvents is approximately 25, 20, and 30 mg/ml, respectively.

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 syringol can be prepared by directly dissolving the solid in aqueous buffers. The solubility of syringol in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Syringol is a phenolic lignan pyrolysis product that has been found in wood smoke.<sup>1,2</sup>

### References

1. Kjällstrand, J. and Petersson, G. Phenolic antioxidants in wood smoke. *Sci. Total Environ.* **277(1-3)**, 69-75 (2001).
2. Wedekind, R., Keski-Rahkonen, P., Robinot, N., et al. Syringol metabolites as new biomarkers for smoked meat intake. *Am. J. Clin. Nutr.* **110(6)**, 1424-1433 (2019).

#### 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, 11/04/2022

#### 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