

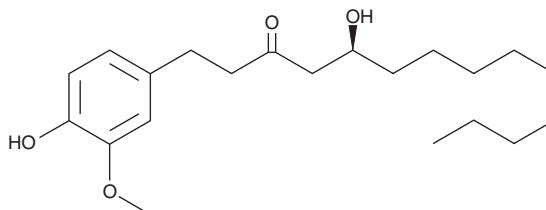
# PRODUCT INFORMATION



## 10-Gingerol

Item No. 11842

CAS Registry No.: 23513-15-7  
Formal Name: 5S-hydroxy-1-(4-hydroxy-3-methoxyphenyl)-3-tetradecanone  
MF: C<sub>21</sub>H<sub>34</sub>O<sub>4</sub>  
FW: 350.5  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 281 nm  
Supplied as: A crystalline solid  
Storage: -20°C  
Stability: ≥4 years  
Item Origin: Plant/*Zingiber officinale*



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

### Laboratory Procedures

10-Gingerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 10-gingerol in the solvent of choice, which should be purged with an inert gas. 10-Gingerol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 10-gingerol in DMSO is approximately 25 mg/ml and approximately 30 mg/ml in ethanol and DMF.

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 10-gingerol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 10-gingerol in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

10-Gingerol is a phenol that has been found in *Z. officinale* and has diverse biological activities.<sup>1-4</sup> It scavenges DPPH (Item No. 14805), superoxide, and hydroxyl radicals in cell-free assays (IC<sub>50</sub>s = 10.47, 1.68, and 1.35 μM, respectively) and inhibits oxidative burst induced by N-formyl-Met-Leu-Phe (fMLP; Item No. 21495) in isolated human polymorphonuclear (PMN) neutrophils when used at a concentration of 6 μM.<sup>1</sup> 10-Gingerol inhibits NETosis and the production of reactive oxygen species (ROS) induced by LPS or lupus-relevant stimuli, including ribonucleoprotein-containing immune complexes and anti-phospholipid antibodies, in isolated human neutrophils.<sup>2</sup> It is active against *M. avium* and *M. tuberculosis* (MICs = 25 and 50 μg/ml, respectively).<sup>3</sup> 10-Gingerol (200 μM) is also cytotoxic to human MDA-MB-231 and MDA-MB-468 and murine 4T1 and EO771 mammary carcinoma cells.<sup>4</sup>

### References

1. Dugasani, S., Pichika, M.R., Nadarajah, V.D., et al. *J. Ethnopharmacol.* **127(2)**, 515-520 (2010).
2. Ali, R.A., Gandhi, A.A., Dai, L., et al. *JCI Insight* **6(3)**, e138385 (2021).
3. Hiserodt, R.D., Franzblau, S.G., and Rosen, R.T. *J. Agric. Food Chem.* **46(7)**, 2504-2508 (1998).
4. Bernard, M.M., McConnery, J.R., and Hoskin, Exp. *Mol. Pathol.* **102(2)**, 370-376 (2017).

#### 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, 12/15/2022

#### 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