

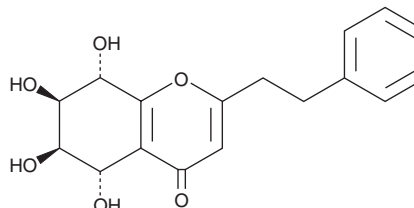
# PRODUCT INFORMATION



## Agarotetrol

Item No. 27662

**CAS Registry No.:** 69809-22-9  
**Formal Name:** (5S,6R,7R,8S)-5,6,7,8-tetrahydro-5,6,7,8-tetrahydroxy-2-(2-phenylethyl)-4H-1-benzopyran-4-one  
**MF:** C<sub>17</sub>H<sub>18</sub>O<sub>6</sub>  
**FW:** 318.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 254 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Lignum Aquilariae Agallochae Resinatum*



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

### Laboratory Procedures

Agarotetrol is supplied as a solid. A stock solution may be made by dissolving the agarotetrol in the solvent of choice, which should be purged with an inert gas. Agarotetrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of agarotetrol in ethanol is approximately 10 mg/ml and approximately 50 mg/ml in DMSO 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 agarotetrol can be prepared by directly dissolving the solid in aqueous buffers. The solubility of agarotetrol in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Agarotetrol is a chromone derivative originally isolated from agarwood, a component of incense and traditional medicines.<sup>1,2</sup>

### References

1. Yoshii, E., Koizumi, T., Oribe, T., *et al.* The structure of agarotetrol, a novel highly oxygenated chromone from agarwood (*jinko*). *Tetrahedron Lett.* **19(41)**, 3921-3924 (1978).
2. Sugiyama, T., Narukawa, Y., Shibata, S., *et al.* Three new 5,6,7,8-tetrahydroxy-5,6,7,8-tetrahydrochromone derivatives enantiomeric to agarotetrol from agarwood. *J. Nat. Med.* **72(3)**, 667-674 (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, 12/19/2022

#### 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