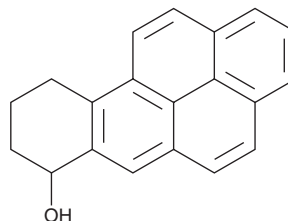


# PRODUCT INFORMATION



## 7,8,9,10-Tetrahydrobenzo[a]pyren-7-ol Item No. 16225

**CAS Registry No.:** 6272-55-5  
**Formal Name:** 7,8,9,10-tetrahydro-benzo[a]pyren-7-ol  
**Synonyms:** 7-HTBP, NSC 30871  
**MF:** C<sub>20</sub>H<sub>16</sub>O  
**FW:** 272.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 246, 267, 278, 314, 328, 345 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

7,8,9,10-Tetrahydrobenzo[a]pyren-7-ol is supplied as a crystalline solid. A stock solution may be made by dissolving the 7,8,9,10-tetrahydrobenzo[a]pyren-7-ol in the solvent of choice, which should be purged with an inert gas. 7,8,9,10-Tetrahydrobenzo[a]pyren-7-ol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 7,8,9,10-tetrahydrobenzo[a]pyren-7-ol in ethanol is approximately 0.2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

7,8,9,10-Tetrahydrobenzo[a]pyren-7-ol is a benzopyrene derivative that is activated by hepatic cytosol into electrophilic sulfuric acid esters, which are capable of forming covalent DNA adducts and inducing mutations.<sup>1,2</sup>

### References

1. Glatt, H., Pauly, K., Frank, H., *et al.* Substance-dependent sex differences in the activation of benzylic alcohols to mutagens by hepatic sulfotransferases of the rat. *Carcinogenesis* **15(11)**, 2605-2611 (1994).
2. Surh, Y.J. and Tannenbaum, S.R. Sulfotransferase-mediated activation of 7,8,9,10-tetrahydro-7-ol, 7,8-dihydrodiol, and 7,8,9,10-tetraol derivatives of benzo[a]pyrene. *Chem. Res. Toxicol.* **8(5)**, 693-698 (1995).

#### 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, 11/17/2022

#### 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