

PRODUCT INFORMATION



Sulisobenzene

Item No. 35225

CAS Registry No.: 4065-45-6

Formal Name: 5-benzoyl-4-hydroxy-2-methoxy-benzenesulfonic acid

Synonyms: Benzophenone 4,
3-Benzoyl-4-hydroxy-6-Methoxybenzenesulfonic Acid,
NSC 60584, UV-284

MF: $C_{14}H_{12}O_6S$

FW: 308.3

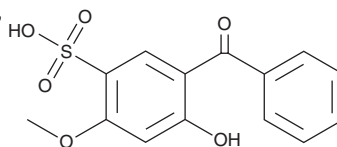
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 225, 243, 287, 325 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Sulisobenzene is supplied as a solid. A stock solution may be made by dissolving the sulisobenzene in the solvent of choice, which should be purged with an inert gas. Sulisobenzene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of sulisobenzene in ethanol is approximately 2 mg/ml and approximately 5 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 sulisobenzene can be prepared by directly dissolving the solid in aqueous buffers. The solubility of sulisobenzene in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Sulisobenzene is a UV filter.¹ Topical application of sulisobenzene reduces UV-induced skin erythema and carcinogenesis in albino mice. Formulations containing sulisobenzene have been used as sunscreens to protect the skin from UVA and UVB rays.

Reference

1. Knox, J.M., Griffin, A.C., and Hakim, R.E. Protection from ultraviolet carcinogenesis. *J. Invest. Dermatol.* **34**, 51-58 (1960).

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, 12/12/2022

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