

# PRODUCT INFORMATION

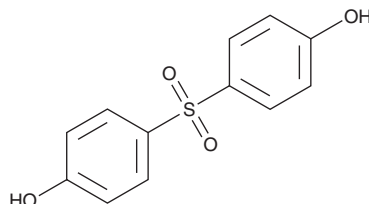


## Bisphenol S

Item No. 45556

**CAS Registry No.:** 80-09-1  
**Formal Name:** 4,4'-sulfonylbis-phenol  
**Synonyms:** Bis(4-hydroxyphenyl)sulfone,  
BPS, NSC 683541, NSC 871,  
4,4'-Sulfonyldiphenol

**MF:** C<sub>12</sub>H<sub>10</sub>O<sub>4</sub>S  
**FW:** 250.3  
**Purity:** ≥98%  
**Supplied as:** A 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

Bisphenol S is supplied as a solid. A stock solution may be made by dissolving the bisphenol S in the solvent of choice, which should be purged with an inert gas. Bisphenol S is soluble (≥10 mg/ml) in ethanol and DMSO.

### Description

Bisphenol S (BPS) is a derivative of bisphenol A (BPA), an environmental contaminant, and an endocrine disruptor.<sup>1,2</sup> It has been detected in surface water and sewage effluent and in personal care products, paper, and food.<sup>3</sup> BPS increases estrogenic activity *in vitro*, in female rats, and in male and female zebrafish. It decreases testosterone levels in male zebrafish and impairs testicular development in male offspring after maternal exposure in rats.<sup>2,3</sup> BPS also has neurotoxic effects, decreasing viability of SK-N-SH cells and primary rat hippocampal neurons, among other cell types.<sup>1</sup> It reduces dendritic length and branching in mice and increases anxiety and impairs learning and memory in rats. BPS increases serum levels of LDL cholesterol and non-esterified fatty acids (NEFA) and induces hepatic lipid accumulation in mice when administered at a dose of 1 mg/kg per day for three months.<sup>4</sup>

### References

1. Cantua, R. and Mulligan, K. Developmental neurotoxicity of bisphenol F and bisphenol S in animal model systems: A literature review. *Neurotoxicology* **108**, 263-280 (2025).
2. Li, B., Tian, Y., Wu, X., *et al.* Maternal bisphenol S exposure impairs testicular development and sperm function in male offspring by disrupting the immune-endocrine network. *Small* e13075 (2026).
3. Rochester, J.R. and Bolden, A.L. Bisphenol S and F: A systematic review and comparison of the hormonal activity of bisphenol A substitutes. *Environ. Health Perspect.* **123**(7), 643-650 (2015).
4. Li, S., Fan, Y., Tang, M., *et al.* Bisphenol S exposure and MASLD: A mechanistic study in mice. *Environ. Health Perspect.* **133**(5), 57009 (2025).

#### 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, 04/23/2026

#### 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