

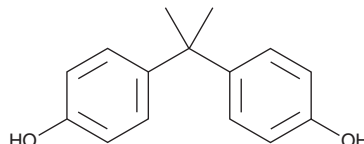
PRODUCT INFORMATION



Bisphenol A

Item No. 36594

CAS Registry No.: 80-05-7
Formal Name: 4,4'-(1-methylethylidene)bis-phenol
Synonyms: BPA, NSC 1767, NSC 17959
MF: $C_{15}H_{16}O_2$
FW: 228.3
Purity: $\geq 98\%$
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

Bisphenol A (BPA) is supplied as a solid. A stock solution may be made by dissolving the BPA in the solvent of choice, which should be purged with an inert gas. BPA is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

BPA is an endocrine disruptor and environmental contaminant.^{1,2} It has estrogenic activity *in vitro*, stimulates uterine, vaginal, and mammary growth and differentiation in rats, and reduces reproduction in zebrafish.^{1,3-5} Prenatal exposure to BPA disrupts metabolic regulation, leading to increased food intake, body and liver weight, and serum insulin levels in adult mice.⁶ BPA has been found in surface water, soil, and air samples, wildlife tissue samples, and human urine.² Formulations containing BPA have been used in the production of plastics, fire retardants, and coatings.

References

1. Matthews, J.B., Twomey, K., and Zacharewski, T.R. In vitro and in vivo interactions of bisphenol A and its metabolite, bisphenol A glucuronide, with estrogen receptors α and β . *Chem. Res. Toxicol.* **14**(2), 149-157 (2001).
2. Corrales, J., Kristofco, L.A., Steele, W.B., et al. Global assessment of bisphenol A in the environment: Review and analysis of its occurrence and bioaccumulation. *Dose Response* **13**(3), 1559325815598308 (2015).
3. Krishnan, A.V., Stathis, P., Permuth, S.F., et al. Bisphenol-A: An estrogenic substance is released from polycarbonate flasks during autoclaving. *Endocrinology* **132**(6), 2279-2286 (1993).
4. Ben-Jonathan, N. and Steinmetz, R. Xenoestrogens: The emerging story of bisphenol A. *Trends Endocrinol. Metab.* **9**(3), 124-128 (1998).
5. Huang, Q., Liu, Y., Chen, Y., et al. New insights into the metabolism and toxicity of bisphenol A on marine fish under long-term exposure. *Environ. Pollut.* **242**(Pt A), 914-921 (2018).
6. Angle, B.M., Do, R.P., Ponzi, D., et al. Metabolic disruption in male mice due to fetal exposure to low but not high doses of bisphenol A (BPA): Evidence for effects on body weight, food intake, adipocytes, leptin, adiponectin, insulin and glucose regulation. *Reprod. Toxicol.* **42**, 256-268 (2013).

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, 07/31/2026

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