

PRODUCT INFORMATION



TBBPA monosulfate

Item No. 9004565

CAS Registry No.: 902525-22-8

Formal Name: 2,6-dibromo-4-[1-(3,5-dibromo-4-hydroxyphenyl)-1-methylethyl]-phenol 1-(hydrogen sulfate)

Synonyms: TBBPA-sulfate, Tetrabromobisphenol A sulfate

MF: $C_{15}H_{12}Br_4O_5S$

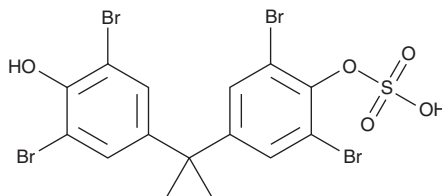
FW: 623.9

Purity: $\geq 98\%$

Supplied as: A neat solid

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

Stability: ≥ 2 years



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

Laboratory Procedures

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

Description

TBBPA monosulfate is an agonist of PPAR γ and an active metabolite of the flame retardant tetrabromobisphenol A ^{1,2} It increases PPAR γ activity in a reporter assay using HGELN cells expressing PPAR γ .

References

1. Riu, A., Maire, A.I., Grimaldi, M., *et al.* Characterization of novel ligands of ER α , ER β , and PPAR γ : The case of halogenated bisphenol A and their conjugated metabolites. *Toxicol. Sci.* **122**(2), 372-382 (2011).
2. Schauer, U.M.D., Völkel, W., and Dekant, W. Toxicokinetics of tetrabromobisphenol a in humans and rats after oral administration. *Toxicol. Sci.* **91**(1), 49-58 (2006).

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, 12/03/2024

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