



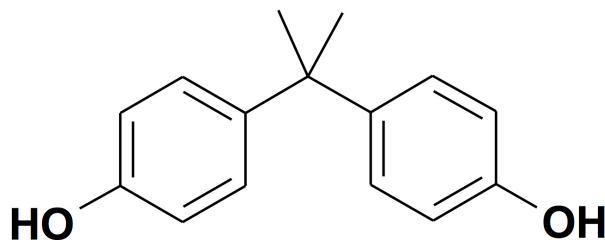
[³H] Bisphenol A

First synthesized in the late 19th century, the worldwide production of bisphenol A (BPA) steadily reached millions of tons by the 1980's. This accelerated manufacturing surge was the direct result of its widespread use in the plastics industry, especially for food and beverage containers. However, at about the same time, growing concern over the safety of BPA also dramatically increased. Mounting evidence pointed to its exposure to and impact on the environment,¹ as well as its multiple routes of human entry through food, inhalation and dermal pathways.² Numerous studies have also documented its role as an endocrine disrupter, acting largely as an estrogenic agent.³ Clearly, BPA is an important and controversial substance with more careful future research required to fully understand it.

To support this critical and emerging research, PerkinElmer now also offers [³H] Bisphenol A (NET1207) at high specific activity and radiopurity. Use of this valuable radioligand will no doubt advance important research in this area.^{4,5,6}

By using radioactive isotopes to directly replace non-radioactive atoms, the biology of the substance you are studying is not altered. The use of radiochemicals is of critical importance in the drug development process for use as radioligands in lead discovery, as metabolic tracers in development, and ADME-Tox studies.

Bisphenol A



References

1. S. Flint, T. Markle, S. Thompson, E. Wallace, J. *Environmental Management*, **2012**, *104*, 19-34.
2. T. Geens, D. Aerts, B. Dominique, C. Berthot, J.-P. Bourguignon, L. Goeyens, P. Lecomte, G. Maghuin-Rogister, A.-M. Pironnet, L. Pussemier, M.-L. Scippo, J. Van Loco, A. Covaci, *Food and Chemical Toxicology*, **2012**, *50*, 3725-3740.
3. P. Alonso-Magdalena, A.B. Ropero, S. Soriano, M. Garcia-Arevalo, C. Ripoll, E. Fuentes, I. Quesada, A. Nadal, *Molecular and Cellular Endocrinology*, **2012**, *355*, 201-207.
4. J.A. Taylor, F.S. vom Saal, W.V. Welshons, B. Drury, G. Rottinghaus, P.A. Hunt, P.-L. Toutain, C.M. Laffont, C.A. Vande Voort, *Environmental Health Perspectives*, **2011**, *119*, 422-430.
5. J.-B. Fini, L. Dolo, J.-P. Cravedi, B. Demeneix, D. Zalko, *Annals of the New York Academy of Sciences*, **2009**, *1163*, 394-397.
6. J.A. Taylor, W.V. Welshons, F.S. vom Saal, *Reproductive Toxicology*, **2008**, *25*, 169-176.