

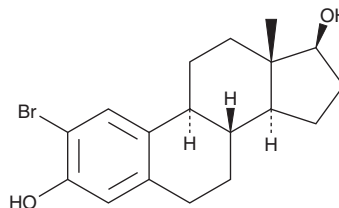
# PRODUCT INFORMATION



## 2-bromo 17 $\beta$ -Estradiol

Item No. 9002145

**CAS Registry No.:** 15833-07-5  
**Formal Name:** 2-bromo-estra-1,3,5(10)-triene-3,17 $\beta$ -diol  
**Synonyms:** 2-bromo E<sub>2</sub>, 2-bromo  $\beta$ -Estradiol, 2-bromo Estradiol, NSC 105460, 2-bromo 17 $\beta$ -Oestradiol  
**MF:** C<sub>18</sub>H<sub>23</sub>BrO<sub>2</sub>  
**FW:** 351.3  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

2-bromo 17 $\beta$ -Estradiol is supplied as a solid. A stock solution may be made by dissolving the 2-bromo 17 $\beta$ -estradiol in the solvent of choice, which should be purged with an inert gas. 2-bromo 17 $\beta$ -Estradiol is soluble in methanol and DMSO.

### Description

2-bromo 17 $\beta$ -Estradiol is a synthetic steroid and an inhibitor of estrogen 2-hydroxylase.<sup>1</sup> It inhibits estrogen 2-hydroxylase in rat liver microsomes when used at concentrations of 50 and 100  $\mu$ M. 2-bromo 17 $\beta$ -Estradiol (5 mg/kg) prevents estrogen-induced inhibition of apomorphine-induced climbing behavior in mice.<sup>2</sup>

### References

1. Brueggemeier, R.W. Inhibition of estrogen 2-hydroxylase. *J. Steroid. Biochem.* **19(5)**, 1683-1685 (1983).
2. Fung, Y.K., Stevens, J.M., Palmer, C.W., Jr., *et al.* Inhibition by bromoestrogens of the effects of estradiol on apomorphine-induced climbing behavior. *Steroids* **49(4-5)**, 287-294 (1987).

#### 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, 11/09/2022

#### 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