PRODUCT INFORMATION



17β-Estradiol-d₂

Item No. 9002846

CAS Registry No.:	53866-33-4	
Formal Name:	estra-1,3,5(10)-triene-2,4-d ₂ -3,17β-diol	
Synonyms:	2,4-Dideuteriostradiol, E_2 - d_2 , Estradiol- d_2 ,	ОН
	β-Estradiol-d ₂ , 17 $β$ -Oestradiol-d ₂ ,	
MF:	$C_{18}H_{22}D_{2}O_{2}^{-1}$	
FW:	274.4	
Chemical Purity:	≥98% (17β-Estradiol)	
Deuterium		
Incorporation:	≥99% deuterated forms (d ₁ -d ₂); ≤1% d ₀	
UV/Vis.:	λ _{max} : 281 nm	
Supplied as:	A crystalline solid	D
Storage:	-20°C	
Stability:	≥4 years	
Information remanant	the preduct exections. Datch execting and the last	ulto are provided on each contificate of analysi

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

Laboratory Procedures

 17β -Estradiol-d₂ contains two deuterium atoms at the 2 and 4 positions. It is intended for use as an internal standard for the quantification of 17β-estradiol (Item Nos. ISO60155 | 10006315) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

17β-Estradiol-d₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the 17β-estradiol-d₂ in the solvent of choice. 17β-Estradiol-d₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 17β -estradiol-d₂ in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

 17β -Estradiol is the major pre-menopausal bioactive estrogen and an active metabolite of testosterone.¹⁻³ It is produced primarily by the ovary, and to a lesser extent, by the adrenals, testes, and adipose tissue, as well as by the placenta during pregnancy.¹ 17β -Estradiol binds to estrogen receptors localized to the nucleus, cytoplasm, and plasma membrane and induces gene transcription or non-genomic intracellular signaling in a location-dependent manner.² It regulates reproductive development in females and spermatogenesis in males, as well as various non-reproductive processes including lipid and glucose homeostasis, bone mass maintenance, cognition, energy balance, and dilation of blood vessels.^{2,3} 17β -Estradiol deficiency due to hypogonadism or 17α -hydroxylase/17,20-lyase deficiency is associated with fatigue, incontinence, osteoporosis, depression, emotional instability, and hot flashes.

References

- 1. Mahboobifard, F., Pourgholami, M.H., Jorjani, M., et al. Estrogen as a key regulator of energy homeostasis and metabolic health. Biomed. Pharmacother. 156, 113808 (2022).
- 2. Kumar, A., Banerjee, A., Sigh, D., et al. Estradiol: A steroid with multiple facets. Horm. Metab. Res. 50(5), 359-374 (2018)
- 3. Sabah-Negah, S., Hajali, V., Moradi, H.R., et al. The impact of estradiol on neurogenesis and cognitive functions in Alzheimer's disease. Cell Mol. Neurobiol. 40(3), 283-299 (2019).

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

uyer 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, 06/22/2023

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