

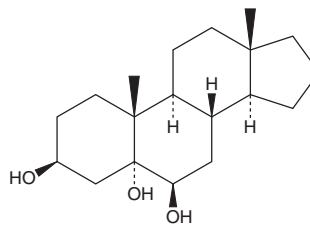
# PRODUCT INFORMATION



## Annaosanchun

Item No. 37572

CAS Registry No.: 4725-51-3  
Formal Name: androstane-3 $\beta$ ,5 $\alpha$ ,6 $\beta$ -triol  
MF: C<sub>19</sub>H<sub>32</sub>O<sub>3</sub>  
FW: 308.5  
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

Annaosanchun is supplied as a solid. A stock solution may be made by dissolving the annaosanchun in the solvent of choice, which should be purged with an inert gas. Annaosanchun is soluble in acetonitrile and chloroform.

### Description

Annaosanchun is a synthetic sterol and neuroprotectant agent.<sup>1</sup> It increases the viability of, and the levels of ATP and the mitochondrial membrane potential in, hypoxia-exposed and reoxygenated primary rat cortical neurons when used at concentrations of 0.1, 0.3, and 1  $\mu$ M. Annaosanchun (10  $\mu$ M) prevents glutamate-induced increases in intracellular acidification, lactate levels, and glycolysis rate, as well as inhibits glutamate-induced apoptosis, in primary rat cerebral granule neurons.<sup>2</sup> It reduces neurological and neuromuscular deficits and infarct volume in a rat model of ischemia-reperfusion injury induced by transient middle cerebral artery occlusion (MCAO) when administered at doses of 0.5, 2, or 10 mg/kg.

### References

1. Chen, J., Leng, T., Chen, W., *et al.* A synthetic steroid 5 $\alpha$ -androst-3 $\beta$ ,5,6 $\beta$ -triol blocks hypoxia/reoxygenation-induced neuronal injuries via protection of mitochondrial function. *Steroids* **78(10)**, 996-1002 (2013).
2. Xue, D., Wei, C., Zhou, Y., *et al.* TRIOL inhibits rapid intracellular acidification and cerebral ischemic injury: The role of glutamate in neuronal metabolic reprogramming. *ACS Chem. Neurosci.* **13(14)**, 2110-2121 (2022).

#### 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, 02/15/2024

#### 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