

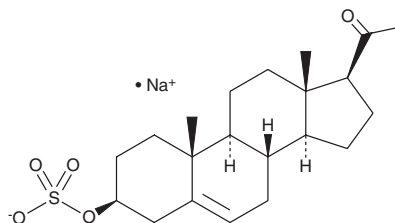
PRODUCT INFORMATION



Pregnenolone sulfate (sodium salt)

Item No. 21004

CAS Registry No.: 1852-38-6
Formal Name: 3-(sulfooxy)-pregn-5-en-20-one, monosodium salt
MF: C₂₁H₃₁O₅S • Na
FW: 418.5
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Pregnenolone sulfate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the pregnenolone sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Pregnenolone sulfate (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pregnenolone sulfate (sodium salt) in ethanol is approximately 2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Pregnenolone sulfate (sodium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pregnenolone sulfate (sodium salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pregnenolone sulfate (sodium salt) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pregnenolone sulfate is a metabolite of the natural steroid hormone pregnenolone (Item No. 19864). Pregnenolone sulfate modulates NMDA receptor responses to exogenously applied glutamate and stimulates transient receptor potential melastatin 3 (TRPM3).¹⁻³

References

1. Sedláček, M., Kořínek, M., Petrovič, M., *et al.* Neurosteroid modulation of ionotropic glutamate receptors and excitatory synaptic transmission. *Physiol. Res.* **57(Suppl 3)**, S49-S57 (2008).
2. Smith, C.C., Gibbs, T.T., and Farb, D.H. Pregnenolone sulfate as a modulator of synaptic plasticity. *Psychopharmacology (Berl)* **231(17)**, 3537-3556 (2014).
3. Thiel, G., Müller, I., and Rössler, O.G. Signal transduction via TRPM3 channels in pancreatic β -cells. *J. Mol. Endocrinol.* **50(3)**, R75-R83 (2013).

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, 12/01/2022

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