

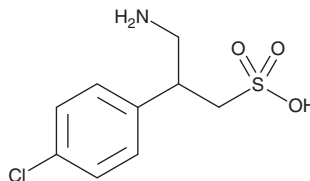
# PRODUCT INFORMATION



## Saclofen

Item No. 17355

**CAS Registry No.:** 125464-42-8  
**Formal Name:**  $\beta$ -(aminomethyl)-4-chloro-benzeethanesulfonic acid  
**MF:** C<sub>9</sub>H<sub>12</sub>ClNO<sub>3</sub>S  
**FW:** 249.7  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 220 nm  
**Supplied as:** A crystalline 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

Saclofen is supplied as a crystalline solid. A stock solution may be made by dissolving the saclofen in the solvent of choice, which should be purged with an inert gas. Saclofen is soluble in DMSO at a concentration of approximately 0.3 mg/ml. It is also soluble at approximately 10 mg/ml in 0.1 M NaOH. We do not recommend storing the aqueous solution for more than one day.

### Description

Saclofen is a sulfonic analog of the inhibitory neurotransmitter  $\gamma$ -aminobutyric acid (GABA) that acts as a competitive antagonist of the GABA<sub>B</sub> receptor ( $IC_{50} = 7.8 \mu\text{M}$ ).<sup>1</sup> This compound can be used to determine the functional roles for the GABA<sub>B</sub> receptor as a mediator of slow inhibitory postsynaptic potentials in the brain.

### Reference

1. Drew, C.A., Johnston, G.A.R., Kerr, D.I.B., *et al.* Inhibition of baclofen binding to rat cerebellar membranes by phaclofen, saclofen, 3-aminopropylphosphonic acid and related GABA<sub>B</sub> receptor antagonists. *Neurosci. Lett.* **113**(1), 107-110 (1990).

#### 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, 10/31/2022

#### 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