

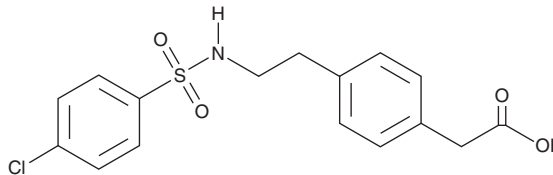
# PRODUCT INFORMATION



## Daltroban

Item No. 14061

**CAS Registry No.:** 79094-20-5  
**Formal Name:** 4-[2-[[[4-chlorophenyl)sulfonyl]amino]ethyl]-benzeneacetic acid  
**Synonyms:** BM 13505, SKF 96148  
**MF:** C<sub>16</sub>H<sub>16</sub>ClNO<sub>4</sub>S  
**FW:** 353.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222 nm  
**Supplied as:** A 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

Daltroban is supplied as a solid. A stock solution may be made by dissolving the daltroban in the solvent of choice, which should be purged with an inert gas. Daltroban is soluble in DMSO.

### Description

Daltroban is a thromboxane receptor (TP) antagonist.<sup>1</sup> It inhibits aggregation of isolated human platelets induced by the TP agonist U-46619 (Item no. 16450; IC<sub>50</sub> = 77 nM). Daltroban increases mean pulmonary arterial pressure (MPAP), mean arterial pressure (MAP), and hematocrit levels in anesthetized rats (ED<sub>50</sub>s = 20, 217, and 94 μg/kg, respectively). It maintains post-reinfusion MAP and superior mesenteric artery flow in a cat model of hemorrhagic shock when administered at a dose of 1 mg/kg.<sup>2</sup> Daltroban (0.25-2 mg/kg) increases survival and decreases circulating platelet count in a rabbit model of sodium arachidonate-induced sudden death.<sup>3</sup>

### References

1. Bertolino, F., Valentin, J.P., Maffre, M., *et al.* Intrinsic activity of the non-prostanoid thromboxane A<sub>2</sub> receptor antagonist, daltroban (BM 13,505), in human platelets *in vitro* and in the rat vasculature *in vivo*. *Br. J. Pharmacol.* **115**(1), 210-216 (1995).
2. Bitterman, H., Yanagisawa, A., and Lefer, A.M. Beneficial actions of thromboxane receptor antagonism in hemorrhagic shock. *Circ. Shock* **20**(1), 1-11 (1986).
3. Lefer, D.J., Mentley, R.K., and Lefer, A.M. Protective effects of a new specific thromboxane antagonist in arachidonate-induced sudden death. *Arch. Int. Pharmacodyn. Ther.* **287**(1), 89-95 (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, 11/02/2022

#### 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