

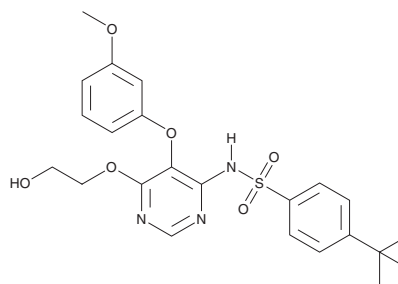
# PRODUCT INFORMATION



Ro 46-2005

Item No. 43523

**CAS Registry No.:** 150725-87-4  
**Formal Name:** 4-(1,1-dimethylethyl)-N-[6-(2-hydroxyethoxy)-5-(3-methoxyphenoxy)-4-pyrimidinyl]-benzenesulfonamide  
**MF:** C<sub>23</sub>H<sub>27</sub>N<sub>3</sub>O<sub>6</sub>S  
**FW:** 473.5  
**Purity:** ≥98%  
**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

Ro 46-2005 is supplied as a solid. A stock solution may be made by dissolving the Ro 46-2005 in the solvent of choice, which should be purged with an inert gas. Ro 46-2005 is soluble (≥10 mg/ml) in DMSO.

## Description

Ro 46-2005 is an antagonist of endothelin receptor type A (ET<sub>A</sub>) and ET<sub>B</sub> (IC<sub>50</sub>s = 220 and 160 nM, respectively).<sup>1</sup> It inhibits endothelin-1-induced secretion of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) in primary rat mesangial cells (IC<sub>50</sub> = 1.8 μM). Ro 46-2005 prevents sarafotoxin S6-induced relaxation in precontracted rat mesenteric arteries (pA<sub>2</sub> = 6.47).<sup>1</sup> It increases cerebral blood flow in a rat model of subarachnoid hemorrhage when administered at a dose of 3 mg/kg. Ro 46-2005 (30 or 100 mg/kg) decreases systolic blood pressure in sodium-depleted monkeys.

## References

1. Breu, V., Löffler, B.-M., and Clozel, M. In vitro characterization of Ro 46-2005, a novel synthetic non-peptide endothelin antagonist of ET<sub>A</sub> and ET<sub>B</sub> receptors. *FEBS Lett.* **334(2)**, 210-214 (1993).
2. Clozel, M., Breu, V., Burri, K., et al. Pathophysiological role of endothelin revealed by the first orally active endothelin receptor antagonist. *Nature* **365(6448)**, 759-761 (1993).

### 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, 06/27/2025

## 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