

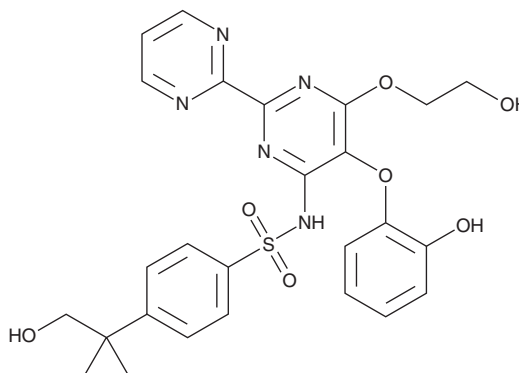
PRODUCT INFORMATION



Hydroxy Desmethyl Bosentan

Item No. 34158

CAS Registry No.: 253688-62-9
Formal Name: 4-(2-hydroxy-1,1-dimethylethyl)-N-[6-(2-hydroxyethoxy)-5-(2-hydroxyphenoxy)[2,2'-bipyrimidin]-4-yl]-benzenesulfonamide
Synonym: Ro 64-1056
MF: C₂₆H₂₇N₅O₇S
FW: 553.6
Purity: ≥90%
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

Hydroxy desmethyl bosentan is supplied as a solid. A stock solution may be made by dissolving the hydroxy desmethyl bosentan in the solvent of choice, which should be purged with an inert gas. Hydroxy desmethyl bosentan is soluble in organic solvents such as methanol and DMSO.

Description

Hydroxy desmethyl bosentan is a metabolite of the endothelin receptor antagonist bosentan (Item No. 11731).¹

Reference

1. van Giersbergen, P.L.M., Gnerre, C., Treiber, A., *et al.* Bosentan, a dual endothelin receptor antagonist, activates the pregnane X nuclear receptor. *Br. J. Clin. Pharmacol.* **54**(5), 561-562 (2002).

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, 11/18/2022

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