

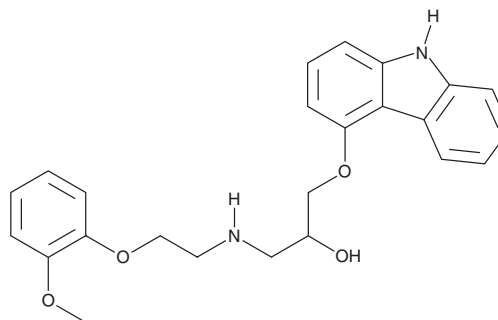
# PRODUCT INFORMATION



## Carvedilol

Item No. 26286

**CAS Registry No.:** 72956-09-3  
**Formal Name:** 1-(9H-carbazol-4-yloxy)-3-[[2-(2-methoxyphenoxy)ethyl]amino]-2-propanol  
**Synonym:** BM 14190  
**MF:** C<sub>24</sub>H<sub>26</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 406.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 244, 286, 320, 333 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Carvedilol (Item No. 26286) is an analytical reference standard categorized as a β-adrenergic receptor antagonist and vasodilator.<sup>1,2</sup> Formulations containing carvedilol have been used to enhance physical performance in athletes.<sup>3</sup> This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 15418).

### References

1. Nichols, A.J., Sulpizio, A.C., Ashton, D.J., *et al.* *In vitro* pharmacologic profile of the novel beta-adrenoceptor antagonist and vasodilator, carvedilol. *Pharmacology* **39(5)**, 327-336 (1989).
2. Bristow, M.R., Roden, R.I., Lowes, B.D., *et al.* The role of third-generation beta-blocking agents in chronic heart failure. *Clin. Cardiol.* **21(12 Suppl 1)**, I3-I13 (1998).
3. Myung, S.-W. and Jo, C.-H. Gas chromatograph-mass spectrometric method for the determination of carvedilol and its metabolites in human urine. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **822(1-2)**, 70-77 (2005).

#### 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, 07/03/2023

#### 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