

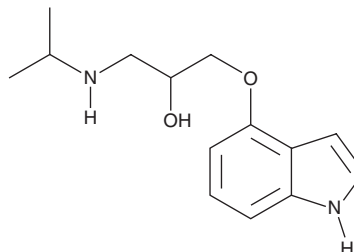
# PRODUCT INFORMATION



## Pindolol

Item No. 21445

<b>CAS Registry No.:</b>	13523-86-9
<b>Formal Name:</b>	1-(1H-indol-4-yloxy)-3-[(1-methylethyl)amino]-2-propanol
<b>Synonym:</b>	DL-Pindolol
<b>MF:</b>	C <sub>14</sub> H <sub>20</sub> N <sub>2</sub> O <sub>2</sub>
<b>FW:</b>	248.3
<b>Purity:</b>	≥99%
<b>UV/Vis.:</b>	λ <sub>max</sub> : 218, 288 nm
<b>Supplied as:</b>	A crystalline solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥2 years



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

### Laboratory Procedures

Pindolol is supplied as a crystalline solid. A stock solution may be made by dissolving the pindolol in the solvent of choice. Pindolol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of pindolol in ethanol is approximately 5 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Pindolol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pindolol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pindolol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pindolol is an antagonist of  $\beta_1$ - and  $\beta_2$ -adrenergic receptors ( $\beta_1$ - and  $\beta_2$ -ARs;  $K_i$ s = 2.6 and 4.8 nM, respectively) and a partial agonist of the  $\beta_3$ -AR, stimulating adenylyl cyclase in membranes of cells expressing the human receptor.<sup>1,2</sup> It is also an antagonist of the serotonin 5-HT<sub>1A</sub> receptor ( $K_i$  = 81.1 nM for inhibition of 5-HT-stimulated GTP $\gamma$ S binding).<sup>3</sup> Pindolol inhibits isoproterenol-induced tachycardia in anesthetized cats ( $ED_{50}$  = 1.8  $\mu$ g/kg).<sup>4</sup> It also decreases mean blood pressure in conscious spontaneously hypertensive rats when administered at a dose of 30 mg/kg per day, but concomitantly increases heart rate.<sup>5</sup> Formulations containing pindolol have been used in the treatment of hypertension.

### References

1. Hicks, P.E., Cavero, I., Manoury, P., et al. *J. Pharmacol. Exp. Ther.* **242**(3), 1025-1034 (1987).
2. Hoffmann, C., Leitz, M.R., Oberdorf-Maass, S., et al. *Naunyn-Schmiedeberg's Arch. Pharmacol.* **369**(2), 151-159 (2004).
3. Krushinski, J.H., Jr., Schaus, J.M., Thompson, D.C., et al. *Bioorg. Med. Chem. Lett.* **17**(20), 5600-5604 (2007).
4. Lubawski, I. and Wale, J. *Eur. J. Pharmacol.* **6**(3), 345-348 (1969).
5. Antonaccio, M.J., High, J., DeForrest, J.M., et al. *J. Pharmacol. Exp. Ther.* **238**(1), 378-387 (1986).

#### 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, 04/18/2023

#### 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