

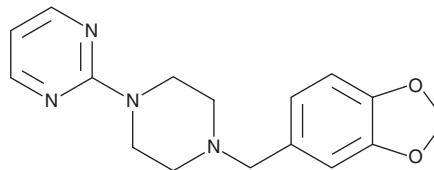
# PRODUCT INFORMATION



## Piribedil

Item No. 30321

**CAS Registry No.:** 3605-01-4  
**Formal Name:** 2-[4-(1,3-benzodioxol-5-ylmethyl)-1-piperazinyl]-pyrimidine  
**MF:** C<sub>16</sub>H<sub>18</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 298.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 242 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

Piribedil is supplied as a solid. A stock solution may be made by dissolving the piribedil in the solvent of choice, which should be purged with an inert gas. Piribedil is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of piribedil in ethanol and DMSO is approximately 5 mg/ml and approximately 10 mg/ml in DMF.

Piribedil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, piribedil should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Piribedil has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Piribedil is a dopamine receptor agonist.<sup>1-4</sup> It binds to dopamine D<sub>2</sub>, D<sub>3</sub>, and D<sub>4</sub> receptors (IC<sub>50</sub>s = 1-1,000 nM in radioligand binding assays).<sup>1</sup> Piribedil (1 and 10 mg/kg) enhances spontaneous object recognition in young adult rats and attenuates age-related memory impairments in the radial arm maze in mice.<sup>2</sup> It reduces immobility time in the forced swim test in mice when administered at doses ranging from 0.63 to 10 mg/kg.<sup>3</sup> Piribedil (0.3 mg/kg) decreases the number of premature and delayed responses in a response time test, indicating improved attentional function, in a rat model of Parkinson's disease induced by 6-OHDA (Item No. 25330) when administered in combination with L-DOPA (Item No. 13248).<sup>4</sup>

### References

1. Cagnotto, A., Parotti, L., and Mennini, T. In vitro affinity of piribedil for dopamine D<sub>3</sub> receptor subtypes, an autoradiographic study. *Eur. J. Pharmacol.* **313**(1-2), 63-67 (1996).
2. Marighetto, A., Valerio, S., Philippin, J.N., et al. Comparative effects of the dopaminergic agonists piribedil and bromocriptine in three different memory paradigms in rodents. *J. Psychopharmacol.* **22**(5), 511-521 (2008).
3. Brocco, M., Dekeyne, A., Papp, M., et al. Antidepressant-like properties of the anti-Parkinson agent, piribedil, in rodents: Mediation by dopamine D<sub>2</sub> receptors. *Behav. Pharmacol.* **17**(7), 559-572 (2006).
4. Turle-Lorenzo, N., Maurin, B., Puma, C., et al. The dopamine agonist piribedil with L-DOPA improves attentional dysfunction: Relevance for Parkinson's disease. *J. Pharmacol. Exp. Ther.* **319**(2), 914-923 (2006).

#### 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, 11/17/2022

#### 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