

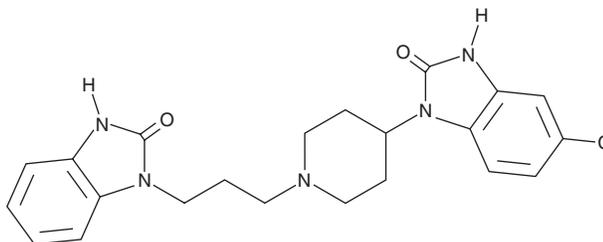
# PRODUCT INFORMATION



## Domperidone

Item No. 18875

**CAS Registry No.:** 57808-66-9  
**Formal Name:** 5-chloro-1-[1-[3-(2,3-dihydro-2-oxo-1H-benzimidazol-1-yl)propyl]-4-piperidinyl]-1,3-dihydro-2H-benzimidazol-2-one  
**Synonyms:** KW 5338, NSC 299589  
**MF:** C<sub>22</sub>H<sub>24</sub>ClN<sub>5</sub>O<sub>2</sub>  
**FW:** 425.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 287 nm  
**Supplied as:** A crystalline 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

Domperidone is supplied as a crystalline solid. A stock solution may be made by dissolving the domperidone in the solvent of choice, which should be purged with an inert gas. Domperidone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of domperidone in these solvents is approximately 10 mg/ml.

Domperidone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, domperidone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Domperidone 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

Domperidone is a dopamine D<sub>2</sub> receptor antagonist (K<sub>i</sub> = 0.3 nM in CHO cells expressing the rat receptor).<sup>1,2</sup> It is selective for dopamine D<sub>2</sub> over D<sub>3</sub> receptors (K<sub>i</sub> = 9.5 nM).<sup>1</sup> Domperidone (0.5-5 μg/kg) inhibits dipropyl dopamine-induced femoral vasodilation in dogs, indicating dopamine D<sub>2</sub> receptor antagonist activity, and has no effect on dopamine-induced vasodilation in the renal vascular bed in dogs when administered at doses up to 5 mg/kg, indicating a lack of activity at dopamine D<sub>1</sub> receptors.<sup>2</sup> Domperidone (0.5 mg/kg) prevents dopamine-induced decreases in gastric antral motility induced by pentagastrin (Item No. 28546) in dogs.<sup>3</sup> It inhibits apomorphine-induced emesis in dogs (ED<sub>50</sub> = 0.031 mg/kg, p.o.).<sup>4</sup> Domperidone also increases serum levels of prolactin in male rats.<sup>5</sup>

### References

1. Sokoloff, P., Giros, B., Martres, M.P., et al. Molecular cloning and characterization of a novel dopamine receptor (D<sub>3</sub>) as a target for neuroleptics. *Nature* **137(6289)**, 146-151 (1990).
2. Kohli, J.D., Glock, D., and Goldberg, L.I. Selective DA<sub>2</sub> versus DA<sub>1</sub> antagonist activity of domperidone in the periphery. *Eur. J. Pharmacol.* **89(1-2)**, 137-141 (1983).
3. Bech, K., Hovendal, C.P., and Andersen, D. Effect of dopamine on pentagastrin-stimulated gastric antral motility in dogs with gastric fistula. *Scand. J. Gastroenterol.* **17(1)**, 103-107 (1982).
4. Niemegeers, C.J.E. Antiemetic specificity of dopamine antagonists. *Psychopharmacology (Berl)*. **78(3)**, 210-213 (1982).
5. Meltzer, H.Y., Simonovic, M., and So, R. Effects of a series of substituted benzamides on rat prolactin secretion and <sup>3</sup>H-spiroperone binding to bovine anterior pituitary membranes. *Life Sci.* **32(25)**, 2877-2886 (1983).

#### 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, 11/28/2022

#### 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