

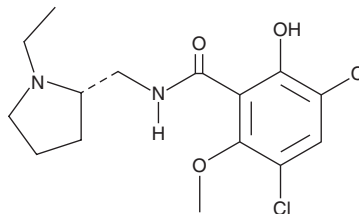
# PRODUCT INFORMATION



## Raclopride

Item No. 17422

**CAS Registry No.:** 84225-95-6  
**Formal Name:** 3,5-dichloro-N-[[[(2S)-1-ethyl-2-pyrrolidinyl]methyl]-2-hydroxy-6-methoxy-benzamide  
**Synonym:** (S)-(-)-Raclopride  
**MF:** C<sub>15</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 347.2  
**Purity:** ≥98%  
**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

Raclopride is supplied as a solid. A stock solution may be made by dissolving the raclopride in the solvent of choice, which should be purged with an inert gas. Raclopride is soluble in organic solvents such as ethanol and DMSO. The solubility of raclopride in these solvents is approximately 20 and 100 mM, respectively.

### Description

Raclopride is a selective dopamine D<sub>2</sub>/D<sub>3</sub> receptor antagonist with K<sub>i</sub> values of 1.8, 3.5, 2,400, and 18,000 nM for D<sub>2</sub>, D<sub>3</sub>, D<sub>4</sub>, and D<sub>1</sub> receptors, respectively.<sup>1</sup> It passes the blood brain barrier and can be used in *in vivo* binding and autoradiography studies of the dopamine system under normal and pathological conditions such as Huntington's disease.<sup>2,3</sup>

### References

1. Seeman, P. and Van Tol, H.H.M. Dopamine receptor pharmacology. *Trends Pharmacol. Sci.* **15**(7), 264-270 (1994).
2. Lidow, M.S., Goldman-Rakic, P.S., Rakic, P., *et al.* Dopamine D<sub>2</sub> receptors in the cerebral cortex: Distribution and pharmacological characterization with [<sup>3</sup>H]raclopride. *Proc. Natl. Acad. Sci. USA* **86**(16), 6412-6416 (1989).
3. Ginovart, N. Imaging the dopamine system with *in vivo* [<sup>11</sup>C]raclopride displacement studies: Understanding the true mechanism. *Mol. Imaging Biol.* **7**(1), 45-52 (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, 01/03/2023

#### 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](http://WWW.CAYMANCHEM.COM)