

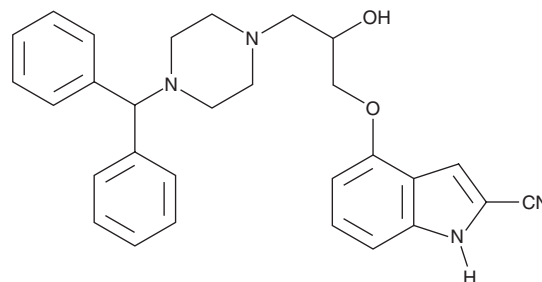
# PRODUCT INFORMATION



## (±)-SDZ-201 106

Item No. 21356

**CAS Registry No.:** 97730-95-5  
**Formal Name:** 4-[3-[4-(diphenylmethyl)-1-piperazinyl]-2-hydroxypropoxy]-1H-indole-2-carbonitrile  
**Synonym:** DPI 201-106  
**MF:** C<sub>29</sub>H<sub>30</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 466.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229, 283 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

(±)-SDZ-201 106 is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-SDZ-201 106 in the solvent of choice, which should be purged with an inert gas. (±)-SDZ-201 106 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)-SDZ-201 106 in ethanol is approximately 1.6 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

(±)-SDZ-201 106 is a diphenylpiperazinylindole derivative that prolongs the open state of cardiac voltage-gated Na<sup>+</sup> channels. It has been shown to increase net Na<sup>+</sup> influx and the activity of reverse mode Na<sup>+</sup>/Ca<sup>2+</sup> exchange, which leads to an increase in intracellular Ca<sup>2+</sup>, as well as to partially inhibit the Na<sup>+</sup> pump in cardiac myocytes.<sup>1</sup> (±)-SDZ-201 106 is also reported to act as an open-channel blocker of delayed-rectifier K<sup>+</sup> channels in NG108-15 neuronal cells.<sup>2</sup>

### References

1. Kaumann, A.H., Richards, D.E., and Russell, D.A. Inhibition of the sodium pump by cardioactive DPI 201-106. *Br. J. Pharmacol.* **91**(1), 3-5 (1987).
2. Wang, Y.-J., Lin, M.-W., Lin, A.-A., *et al.* Evidence for state-dependent block of DPI 201-106, a synthetic inhibitor of Na<sup>+</sup> channel inactivation, on delayed-rectifier K<sup>+</sup> current in pituitary tumor (GH3) cells. *J. Physiol. Pharmacol.* **59**(3), 409-423 (2008).

#### 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, 10/19/2022

#### 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