

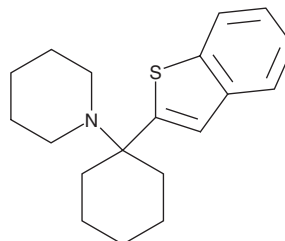
# PRODUCT INFORMATION



## Benocyclidine

Item No. 11737

**CAS Registry No.:** 112726-66-6  
**Formal Name:** 1-(1-benzo[b]thien-2-ylcyclohexyl)-piperidine  
**Synonyms:** BCP, Benzothiophenylcyclohexylpiperidine, BTCP, GK 13  
**MF:** C<sub>19</sub>H<sub>25</sub>NS  
**FW:** 299.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229, 270, 300 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Benocyclidine (BCP) is a derivative of phencyclidine with a benzothiophenyl group instead of a phenyl ring. It acts as a potent and selective dopamine reuptake inhibitor ( $IC_{50} = 8$  nM) with negligible affinity for the NMDA receptor-linked phencyclidine receptor ( $K_{0.5} = 6$  μM).<sup>1,2</sup> BCP has been used to label the dopamine transporter in the mouse brain.<sup>3</sup> This product is intended for forensic and research applications.

### References

1. Vignon, J., Pinet, V., Cerruti, C., *et al.* [<sup>3</sup>H]N-[1-(2-benzo(b)thiophenyl)cyclohexyl]piperidine ([<sup>3</sup>H]BTCP): A new phencyclidine analog selective for the dopamine uptake complex. *Eur. J. Pharmacol.* **148(3)**, 427-436 (1988).
2. Chaudieu, I., Vignon, J., Chicheportiche, M., *et al.* Role of the aromatic group in the inhibition of phencyclidine binding and dopamine uptake by PCP analogs. *Pharmacol. Biochem. Behav.* **32(3)**, 699-705 (1989).
3. Maurice, T., Vignon, J., Kamenka, J.-M., *et al.* *In vivo* labelling of the mouse dopamine uptake complex with the phencyclidine derivative [<sup>3</sup>H]BTCP. *Neurosci. Lett.* **101(2)**, 234-238 (1989).

#### 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, 07/18/2023

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