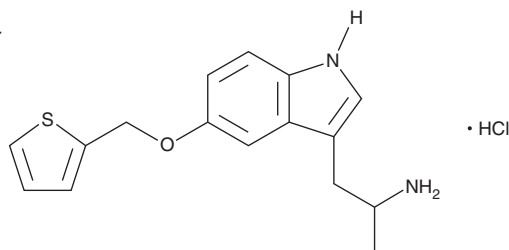


# PRODUCT INFORMATION



**BW 723C86**  
Item No. 70090

**CAS Registry No.:** 160521-72-2  
**Formal Name:**  $\alpha$ -methyl-5-(2-thienylmethoxy)-1H-indole-3-ethanamine, monohydrochloride  
**MF:** C<sub>16</sub>H<sub>18</sub>N<sub>2</sub>OS • HCl  
**FW:** 322.9  
**Purity:** ≥98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 224 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

BW 723C86 is supplied as a crystalline solid. A stock solution may be made by dissolving the BW 723C86 in the solvent of choice, which should be purged with an inert gas. BW 723C86 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BW 723C86 in these solvents is approximately 6, 52, and 55 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of BW 723C86 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of BW 723C86 in PBS (pH 7.2) is approximately 0.16 mg/ml. We do not recommend storing the aqueous solution for more than one day.

## Description

The transduction of neurobehavioral effects by serotonin (5-hydroxy tryptamine; 5-HT) is mediated *via* at least 6 major 5-HT receptor subtypes. Non-selective 5-HT receptor agonists, such as mCPP, induce many behavioral responses in rats, including increased anxiety, hypophagia, and hypoactivity. BW 723C86 is in fact anxiolytic, and its only other prominent behavioral effect is hyperphagia.<sup>1</sup> This effect is likely due to the 10-fold selectivity of BW 723C86 for the HT<sub>2B</sub> receptor.<sup>2</sup> Direct infusion of 1-3  $\mu$ g of BW 723C86 into the cerebral ventricles of rats evokes maximal behavioral responses; the response was attenuated at both higher and lower doses.<sup>2</sup>

## References

1. Kennett, G.A., Bright, F., Trail, B., *et al.* Effects of the 5-HT<sub>2B</sub> receptor agonist, BW 723C86, on three rat models of anxiety. *Br. J. Pharmacol.* **117**(7), 1443-1448 (1996).
2. Kennett, G.A., Ainsworth, K., Trail, B., *et al.* BW 723C86, a 5-HT<sub>2B</sub> receptor agonist, causes hyperphagia and reduced grooming in rats. *Neuropharmacology* **36**(2), 233-239 (1997).

### 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