

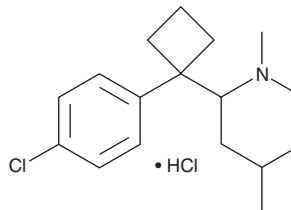
PRODUCT INFORMATION



Sibutramine (hydrochloride)

Item No. 14476

CAS Registry No.: 84485-00-7
Formal Name: 1-(4-chlorophenyl)-N,N-dimethyl- α -(2-methylpropyl)-cyclobutanemethanamine, monohydrochloride
Synonym: BTS-54524
MF: C₁₇H₂₆NCl • HCl
FW: 316.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 224 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Sibutramine (hydrochloride) (Item No. 14476) is an analytical reference standard categorized as a phenylethylamine. It inhibits reuptake of serotonin (Item No. 14332) and norepinephrine (Item No. 16673) via the respective transporters.¹ Sibutramine can be detected in plasma and urine.^{2,3} Formulations containing sibutramine were previously used as anorexiant for obesity but its use was terminated due to adverse cardiovascular events.¹ Sibutramine is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

References

1. Kang, J.G. and Park, C.-Y. Anti-obesity drugs: A review about their effects and safety. *Diabetes Metab. J.* **36(1)**, 13-25 (2012).
2. Jain, D.S., Subbaiah, G., Sanyal, M., et al. Liquid chromatography/electrospray ionization tandem mass spectrometry validated method for the simultaneous quantification of sibutramine and its primary and secondary amine metabolites in human plasma and its application to a bioequivalence study. *Rapid Commun. Mass Spectrom.* **20(23)**, 3509-3521 (2006).
3. Crevelin, E.J., Salami, F.H., Alves, M.N., et al. Direct analysis of amphetamine stimulants in a whole urine sample by atmospheric solids analysis probe tandem mass spectrometry. *J. Am. Soc. Mass Spectrom.* **27(5)**, 944-947 (2016).

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, 06/16/2023

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