

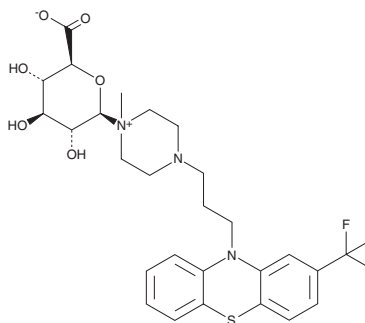
PRODUCT INFORMATION



Trifluoperazine N-Glucuronide

Item No. 34299

CAS Registry No.: 165602-90-4
Formal Name: 1-β-D-glucopyranuronosyl-1-methyl-4-[3-[2-(trifluoromethyl)-10H-phenothiazin-10-yl]propyl]-piperazinium, inner salt
Synonym: TFP N-Glucuronide
MF: C₂₇H₃₂F₃N₃O₆S
FW: 583.6
Purity: ≥70%
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

Trifluoperazine N-glucuronide is supplied as a solid. A stock solution may be made by dissolving the trifluoperazine N-glucuronide in the solvent of choice. Trifluoperazine N-glucuronide is soluble in organic solvent methanol, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Trifluoperazine N-glucuronide is a metabolite of the antipsychotic agent trifluoperazine (Item No. 15068).^{1,2} It is formed from trifluoperazine by the UDP-glucuronosyltransferase (UGT) isoform UGT1A4.²

References

1. Luo, H., Hawes, E.M., McKay, G., *et al.* N⁺-glucuronidation of aliphatic tertiary amines in human: antidepressant versus antipsychotic drugs. *Xenobiotica* **25**(3), 291-301 (1995).
2. Fujiwara, R., Nakajima, M., Yamanaka, H., *et al.* Interactions between human UGT1A1, UGT1A4, and UGT1A6 affect their enzymatic activities. *Drug Metab. Dispos.* **35**(10), 1781-1787 (2007).

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, 11/11/2022

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