

PRODUCT INFORMATION

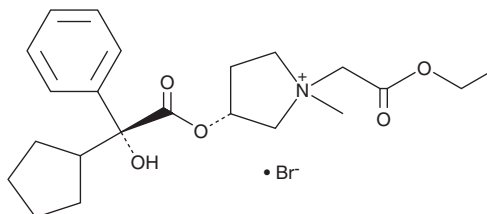


Sofpironium (bromide)

Item No. 38794

CAS Registry No.: 1628106-94-4
Formal Name: (3R)-3-[[[(2R)-2-cyclopentyl-2-hydroxy-2-phenylacetyl]oxy]-1-(2-ethoxy-2-oxoethyl)-1-methylpyrrolidinium, monobromide

Synonym: BBI 4000
MF: C₂₂H₃₂NO₅ • Br
FW: 470.4
Purity: ≥95%
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

Sofpironium (bromide) is supplied as a solid. A stock solution may be made by dissolving the sofpiroonium (bromide) in the solvent of choice, which should be purged with an inert gas. Sofpironium (bromide) is slightly soluble in acetonitrile.

Sofpironium (bromide) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Sofpironium is a derivative of the muscarinic acetylcholine receptor antagonist glycopyrrolate (Item No. 16498).¹ Formulations containing sofpiroonium have been used in the treatment of hyperhidrosis.

Reference

1. Tóth-Sarudy, E., Tóth, G., Pallagi, I., *et al.* Preparation and biological effects of pure stereoisomeric novel soft anticholinergics. *Pharmazie* **61**(2), 90-96 (2006).

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, 12/06/2023

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