

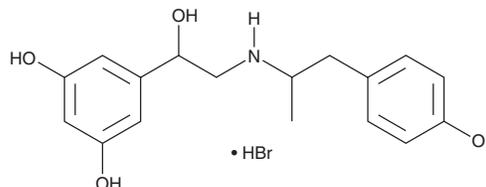
PRODUCT INFORMATION



Fenoterol (hydrobromide)

Item No. 21293

CAS Registry No.: 1944-12-3
Formal Name: 5-[1-hydroxy-2-[[2-(4-hydroxyphenyl)-1-methylethyl]amino]ethyl]-1,3-benzenediol, monohydrobromide
Synonym: Phenoterol
MF: C₁₇H₂₁NO₄ • HBr
FW: 384.3
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 278 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

Fenoterol (Item No. 21293) is an analytical reference standard categorized as a β₂-adrenergic receptor agonist.¹ β₂-adrenergic receptor agonists, including fenoterol, have been used as performance-enhancing drugs in sports doping and are banned by the World Anti-Doping Agency (WADA).¹ This product is intended for research and forensic applications.

Reference

1. Kolmonen, M., Leinonen, A., Pelander, A., *et al.* A general screening method for doping agents in human urine by solid phase extraction and liquid chromatography/time-of-flight mass spectrometry. *Anal. Chim. Acta* **585(1)**, 94-102 (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, 07/06/2023

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