

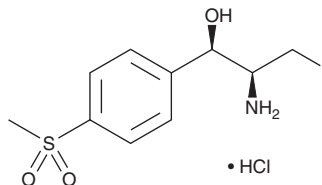
PRODUCT INFORMATION



Florfenicol amine (hydrochloride)

Item No. 18631

CAS Registry No.: 108656-33-3
Formal Name: αR-[(1S)-1-amino-2-fluoroethyl]-4-(methylsulfonyl)-benzenemethanol, monohydrochloride
MF: C₁₀H₁₄FNO₃S • HCl
FW: 283.7
Purity: ≥95%
UV/Vis.: λ_{max}: 222, 265, 272 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

Florfenicol amine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the florfenicol amine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Florfenicol amine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of florfenicol amine (hydrochloride) in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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 florfenicol amine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of florfenicol amine (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Florfenicol amine is a metabolite of the broad-spectrum antibiotic florfenicol (Item No. 19458).¹

Reference

1. Anadón, A., Martínez, M.A., Martínez, M., *et al.* Plasma and tissue depletion of florfenicol and florfenicol-amine in chickens. *J. Agric. Food Chem.* **56**(22), 11049-11056 (2008).

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, 06/20/2023

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