

PRODUCT INFORMATION

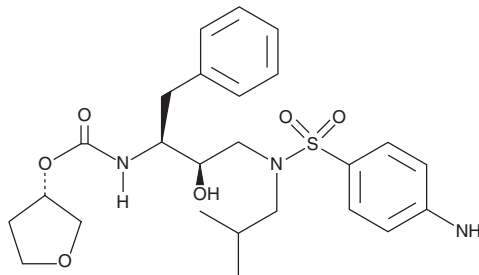


Amprenavir

Item No. 15369

CAS Registry No.: 161814-49-9
Formal Name: N-[(1S,2R)-3-[[[4-aminophenyl]sulfonyl](2-methylpropyl)amino]-2-hydroxy-1-(phenylmethyl)propyl]-carbamic acid, (3S)-tetrahydro-3-furanyl ester

Synonym: VX-478
MF: C₂₅H₃₅N₃O₆S
FW: 505.6
Purity: ≥95%
UV/Vis.: λ_{max}: 268 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

Amprenavir is supplied as a crystalline solid. A stock solution may be made by dissolving the amprenavir in the solvent of choice, which should be purged with an inert gas. Amprenavir is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of amprenavir in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

Amprenavir is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, amprenavir should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Amprenavir has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Amprenavir is an inhibitor of HIV protease ($K_i = 0.04$ nM).¹ It inhibits the cytopathic effects of HIV-1 in MT-4 cells ($IC_{50} = 150$ nM). Formulations containing amprenavir have been used in combination with other antiretroviral agents in the treatment of HIV-1 infection.

Reference

- Sherrill, R.G., Furfine, E.S., Hazen, R.J., *et al.* Synthesis and antiviral activities of novel N-alkoxy-arylsulfonamide-based HIV protease inhibitors. *Bioorg. Med. Chem. Lett.* **15(15)**, 3560-3564 (2005).

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, 11/30/2022

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