

PRODUCT INFORMATION



BAY-1895344 (hydrochloride)

Item No. 26536

Formal Name: 2-[(3R)-3-methyl-4-morpholinyl]-4-(1-methyl-1H-pyrazol-5-yl)-8-(1H-pyrazol-3-yl)-1,7-naphthyridine, monohydrochloride

MF: C₂₀H₂₁N₇O • HCl

FW: 411.9

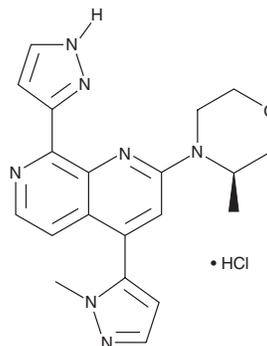
Purity: ≥98%

UV/Vis.: λ_{max}: 256, 296, 431 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

BAY-1895344 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY-1895344 (hydrochloride) in the solvent of choice. BAY-1895344 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. It is also soluble in water. The solubility of BAY-1895344 (hydrochloride) in ethanol, DMSO, and water is approximately 30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BAY-1895344 is an inhibitor of ataxia-telangiectasia and Rad3-related protein (ATR) kinase that inhibits proliferation in a panel of human tumor cell lines (mean IC₅₀ = 78 nM).¹ It halts tumor growth in ovarian and colorectal mouse xenograft models and induces complete tumor remission in mantle cell lymphoma mouse xenograft models. BAY-1895344 also exhibits synergistic anti-tumor activity with radium-223 in a mouse xenograft model of castration-resistant prostate cancer bone metastases.

Reference

1. Wengner, A.M., Siemeister, G., Luecking, U., *et al.* Abstract 836: ATR inhibitor BAY 1895344 shows potent anti-tumor efficacy in monotherapy and strong combination potential with the targeted alpha therapy Radium-223 dichloride in preclinical tumor models. *Cancer Res.* **77(13 Suppl)**, 836-836 (2017).

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, 10/07/2022

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