

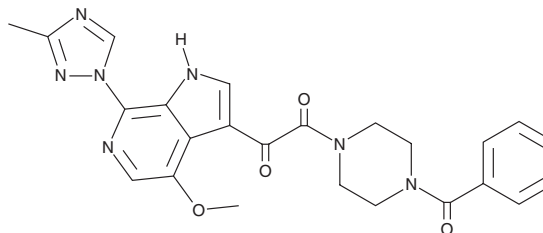
PRODUCT INFORMATION



BMS 626529

Item No. 18413

CAS Registry No.: 701213-36-7
Formal Name: 1-(4-benzoyl-1-piperazinyl)-2-[4-methoxy-7-(3-methyl-1H-1,2,4-triazol-1-yl)-1H-pyrrolo[2,3-c]pyridin-3-yl]-1,2-ethanedione
MF: C₂₄H₂₃N₇O₄
FW: 473.5
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 315 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

BMS 626529 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 626529 in the solvent of choice. BMS 626529 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of BMS 626529 in these solvents is approximately 0.2, 0.3, and 0.1 mg/ml, respectively.

BMS 626529 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMS 626529 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BMS 626529 has a solubility of approximately 0.03 mg/ml in a 1:30 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BMS 626529 is an inhibitor of HIV-1 attachment.^{1,2} It binds to non-ligand bound HIV-1 gp120 to inhibit HIV-1 interaction with host CD4⁺ T cells and subsequent HIV-1 binding and cell entry.¹ BMS 626529 reduces infectivity of laboratory strains and clinical isolates of HIV-1 (EC₅₀s = 0.4-2,000 and 25-2,000 nM, respectively) with cytotoxic concentration (CC₅₀) values greater than 100 μM in a panel of mammalian cell lines.²

References

- Langley, D.R., Kimura, S.R., Sivaprakasam, P., *et al.* Homology models of the HIV-1 attachment inhibitor BMS-626529 bound to gp120 suggest a unique mechanism of action. *Proteins* **83**(2), 331-350 (2015).
- Nowicka-Sans, B., Gong, Y.-F., McAuliffe, B., *et al.* *In vitro* antiviral characteristics of HIV-1 attachment inhibitor BMS-626529, the active component of the prodrug BMS-663068. *Antimicrob. Agents Chemother.* **56**(7), 3498-3507 (2012).

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, 12/02/2022

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