

# PRODUCT INFORMATION

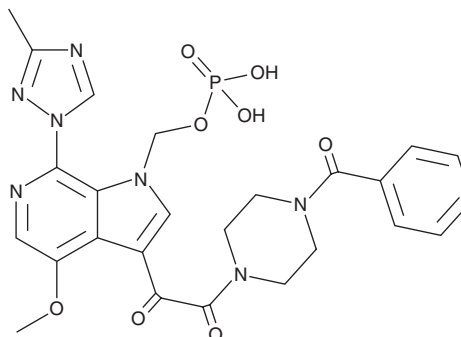


## BMS 663068

Item No. 37708

**CAS Registry No.:** 864953-29-7  
**Formal Name:** 1-(4-benzoyl-1-piperazinyl)-2-[4-methoxy-7-(3-methyl-1H-1,2,4-triazol-1-yl)-1-[(phosphonoxy)methyl]-1H-pyrrolo[2,3-c]pyridin-3-yl]-1,2-ethanedione

**Synonym:** Fostemsavir  
**MF:** C<sub>25</sub>H<sub>26</sub>N<sub>7</sub>O<sub>8</sub>P  
**FW:** 583.5  
**Purity:** ≥98%  
**Supplied as:** A 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 663068 is supplied as a solid. A stock solution may be made by dissolving the BMS 663068 in the solvent of choice, which should be purged with an inert gas. BMS 663068 is soluble in organic solvents such as methanol and DMSO. BMS 663068 is slightly soluble in water and acetonitrile.

BMS 663068 is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

BMS 663068 is a prodrug form of the HIV-1 attachment inhibitor BMS 626529 (Item No. 18413).<sup>1</sup> Formulations containing BMS 663068 have been used in the treatment of multidrug-resistant HIV-1 infections.

### Reference

1. 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, 11/01/2023

#### 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