

# PRODUCT INFORMATION

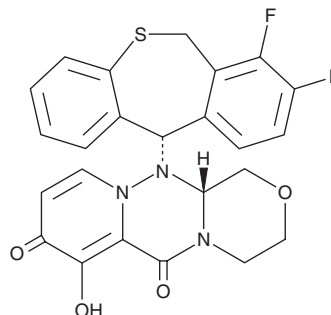


## Baloxavir

Item No. 35125

**CAS Registry No.:** 1985605-59-1  
**Formal Name:** (12aR)-12-[(11S)-7,8-difluoro-6,11-dihydrodibenzo[b,e]thiepin-11-yl]-3,4,12,12a-tetrahydro-7-hydroxy-1H-[1,4]oxazino[3,4-c]pyrido[2,1-f][1,2,4]triazine-6,8-dione

**Synonyms:** BXA, S-033447  
**MF:** C<sub>24</sub>H<sub>19</sub>F<sub>2</sub>N<sub>3</sub>O<sub>4</sub>S  
**FW:** 483.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 260 nm  
**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

Baloxavir is supplied as a solid. A stock solution may be made by dissolving the baloxavir in the solvent of choice, which should be purged with an inert gas. Baloxavir is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of baloxavir in these solvents is approximately 1 and 10 mg/ml, respectively.

### Description

Baloxavir is an active metabolite of baloxavir marboxil (Item No. 33214).<sup>1</sup> It inhibits influenza virus cap-dependent endonuclease (CEN) and CEN/RNA-dependent RNA polymerase (CEN/RdRp) activities in cell-free assays (IC<sub>50</sub>s = 2.5 and 1.6 nM, respectively). Baloxavir reduces viral titers in MDCK cells infected with laboratory, clinical, or zoonotic strains of influenza A or B (EC<sub>50</sub>s = 0.46-6.5 nM).

### Reference

1. Noshi, T., Kitano, M., Taniguchi, K., *et al.* In vitro characterization of baloxavir acid, a first-in-class cap-dependent endonuclease inhibitor of the influenza virus polymerase PA subunit. *Antiviral Res.* **160**, 109-117 (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, 11/14/2022

#### 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