

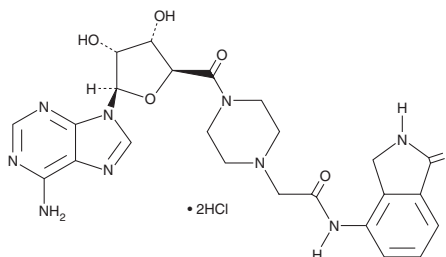
PRODUCT INFORMATION



EB 47 (hydrochloride)

Item No. 34684

CAS Registry No.: 1190332-25-2
Formal Name: 5'-deoxy-5'-[4-[2-[(2,3-dihydro-1-oxo-1H-isoindol-4-yl)amino]-2-oxoethyl]-1-piperazinyl]-5'-oxo-adenosine, dihydrochloride
MF: C₂₄H₂₇N₉O₆ • 2HCl
FW: 610.5
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

EB 47 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the EB 47 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. EB 47 (hydrochloride) is soluble in the organic solvent DMSO (sonicated) at a concentration of approximately 50 mg/ml.

Description

EB 47 is an inhibitor of poly(ADP-ribose) polymerase 1 (PARP1; IC₅₀ = 45 nM) and a dual site inhibitor of tankyrase 2 (TNKS2; IC₅₀ = 45 nM).^{1,2} It also inhibits tankyrase 1 (TNKS1) and PARP10 (IC₅₀s = 410 and 1,179 nM, respectively).^{2,3} EB 47 (10 mg/kg per hour) reduces infarct volume in a rat model of ischemia-reperfusion injury induced by middle cerebral artery occlusion (MCAO).⁴

References

1. Jagtap, P.G., Southan, G.J., Baloglu, E., *et al.* The discovery and synthesis of novel adenosine substituted 2,3-dihydro-1H-isoindol-1-ones: Potent inhibitors of poly(ADP-ribose) polymerase-1 (PARP-1). *Bioorg. Med. Chem. Lett.* **14(1)**, 81-85 (2004).
2. Haikarainen, T., Narwal, M., Joensuu, P., *et al.* Evaluation and structural basis for the inhibition of tankyrases by PARP inhibitors. *ACS Med. Chem. Lett.* **5(1)**, 18-22 (2013).
3. Venkannagari, H., Fallarero, A., Feijs, K.L.H., *et al.* Activity-based assay for human mono-ADP-ribosyltransferases ARTD7/PARP15 and ARTD10/PARP10 aimed at screening and profiling inhibitors. *Eur. J. Pharm. Sci.* **49(2)**, 148-156 (2013).
4. Ferraris, D.V. Evolution of poly(ADP-ribose) polymerase-1 (PARP-1) inhibitors. From concept to clinic. *J. Med. Chem.* **53(12)**, 4561-4584 (2010).

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