

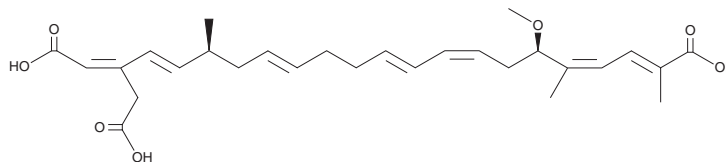
PRODUCT INFORMATION



Bongkreikic Acid

Item No. 38230

CAS Registry No.: 11076-19-0
Formal Name: 20-(carboxymethyl)-6R-methoxy-2,5,17S-trimethyl-2E,4Z,8Z,10E,14E,18E,20Z-docosaheptaenedioic acid
Synonym: Bongkreik Acid
MF: C₂₈H₃₈O₇
FW: 486.6
Purity: ≥98%
Supplied as: A solution in 0.01M Tris buffer pH 7.5
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Pseudomonas gladioli*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Bongkreikic acid is supplied as a solution in 0.01M Tris buffer pH 7.5. To change the solvent, simply evaporate the 0.01M Tris buffer pH 7.5 under a gentle stream of nitrogen and immediately add the solvent of choice. Bongkreikic acid is miscible in DMSO.

Description

Bongkreikic acid is polyketide and bacterial toxin that has been found in *B. gladioli*.^{1,2} It is an inhibitor of ADP/ATP translocase (AAT), a carrier protein that imports ADP into the mitochondrial matrix and exports ATP into the intermembrane space.^{1,3} It prevents ADP import and inhibits mitochondrial respiration ($K_i = 20$ nM in isolated rat liver mitochondria).^{4,5} Bongkreikic acid induces cell death in MCF-7 human breast cancer cells deprived of estradiol for six months ($EC_{50} = 2.58$ μ M).⁶ It is toxic to mice ($LD_{50} = 0.68$ - 6.84 mg/kg, p.o.).⁷

References

1. Henderson, P.J.F. and Lardy, H.A. Bongkreikic acid. An inhibitor of the adenine nucleotide translocase of mitochondria. *J. Biol. Chem.* **245**(6), 1319-1326 (1970).
2. Xie, X., Khosla, C., and Cane, D.E. Elucidation of the stereospecificity of C-methyltransferases from *trans*-at polyketide synthases. *J. Am. Chem. Soc.* **139**(17), 6102-6105 (2017).
3. Lauquin, G.J., Dulpaa, A.M., Klein, G., *et al.* Isobongkreikic acid, a new inhibitor of mitochondrial ADP-ATP transport: Radioactive labeling and chemical and biological properties. *Biochemistry* **15**(11), 2323-2327 (1976).
4. Kligenberg, M., Grebe, K., and Heldt, H.W. On the inhibition of the adenine nucleotide translocation by bongkreikic acid. *Biochem. Biophys. Res. Commun.* **39**(3), 344-351 (1970).
5. Ruprecht, J.J., King, M.S., Zögg, T., *et al.* The molecular mechanism of transport by the mitochondrial ADP/ATP carrier. *Cell* **176**(3), 435-447 (2019).
6. Takeda, S., Okazaki, H., Kudo, T., *et al.* Bongkreikic acid as a Warburg effect modulator in long-term estradiol-deprived MCF-7 breast cancer cells. *Anticancer Res.* **36**(10), 5171-5182 (2016).
7. Hu, W.J., Zhang, G.S., Chu, F.S., *et al.* Purification and partial characterization of flavotoxin A. *Appl. Environ. Microbiol.* **48**(4), 690-693 (1984).

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, 01/16/2024

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