

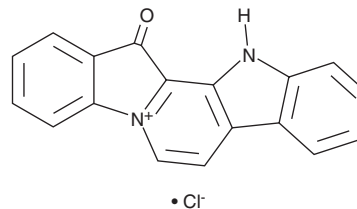
PRODUCT INFORMATION



Fascaplysin (chloride)

Item No. 21715

CAS Registry No.: 114719-57-2
Formal Name: 12,13-dihydro-13-oxo-pyrido[1,2-a:3,4-b']diindol-5-ium, monochloride
Synonym: NSC 622398
MF: C₁₈H₁₁N₂O • Cl
FW: 306.8
Purity: ≥95%
UV/Vis.: λ_{max}: 215, 262, 299, 407 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

Fascaplysin (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the fascaplysin (chloride) in the solvent of choice, which should be purged with an inert gas. Fascaplysin (chloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of fascaplysin (chloride) in these solvents is approximately 16 and 0.14 mg/ml, respectively.

Fascaplysin (chloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fascaplysin (chloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fascaplysin (chloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fascaplysin is an inhibitor of cyclin D kinase 4/ cyclin D1 (IC₅₀ = 0.35 μM) that was originally isolated from the marine sponge *Thorectandra*.¹ It is significantly less selective for Cdk6/cyclin D1 (IC₅₀ = 3.4 μM) and does not inhibit other Cdks and tyrosine kinases.¹ Fascaplysin is reported to have antiproliferative activity against retinoblastoma-positive tumor cells, blocking retinoblastoma phosphorylation and inducing G₁ arrest.¹ Inhibition of Cdk4 *via* fascaplysin has been shown to induce peroxisome-proliferator-activated receptor γ-1α deacetylation (IC₅₀ = 0.7 μM) and has been used to demonstrate a role for insulin-activated cyclinD1-Cdk4 signaling in the control of glucose metabolism.²

References

1. Soni, R., Muller, L., Furet, P., *et al.* Inhibition of cyclin-dependent kinase 4 (Cdk4) by fascaplysin, a marine natural product. *Biochem. Biophys. Res. Commun.* **275**(3), 877-884 (2000).
2. Lee, Y., Dominy, J.E., Choi, Y.J., *et al.* Cyclin D1-Cdk4 controls glucose metabolism independently of cell cycle progression. *Nature* **510**(7506), 547-551 (2014).

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