

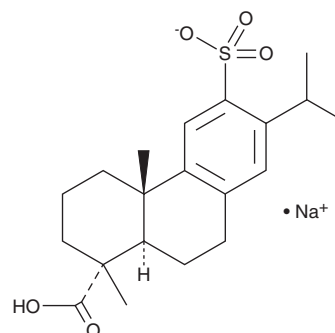
PRODUCT INFORMATION



Ecabet (sodium salt)

Item No. 28013

CAS Registry No.: 86408-72-2
Formal Name: (1R,4aS,10aR)-1,2,3,4,4a,9,10,10a-octahydro-1,4a-dimethyl-7-(1-methylethyl)-6-sulfo-1-phenanthrenecarboxylic acid, monosodium salt
Synonym: TA-2711
MF: C₂₀H₂₇O₅S • Na
FW: 402.5
Purity: ≥98%
UV/Vis.: λ_{max}: 222 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

Ecabet (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the ecabet (sodium salt) in the solvent of choice, which should be purged with an inert gas. Ecabet (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ecabet (sodium salt) in these solvents is approximately 3, 30, and 25 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ecabet (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ecabet (sodium salt) in PBS, pH 7.2, is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ecabet is an antiulcerative and gastroprotective agent.¹⁻⁴ It reduces the viability of *H. pylori* grown under urea-supplemented acidic conditions in a concentration-dependent manner.¹ Ecabet inhibits *H. pylori* urease activity in a pH-dependent manner (IC₅₀s = 2.1-2.6 and >16 mg/ml at pH values of 5 and 8, respectively). It inhibits adhesion of *H. pylori* to MKN-28 cells when used at concentrations of 1 and 2 mg/ml.³ *In vivo*, ecabet (25 and 100 mg/kg, p.o.) increases levels of prostaglandin E₂ (PGE₂; Item No. 14010) in rat gastric mucosa.³ It inhibits formation of hemorrhagic lesions in esophageal mucosa and reduces gastric juice pepsin activity in a rat model of reflux induced by fore-stomach and pyloric ligation when administered at a dose of 30 mg/kg.⁴

References

1. Ito, Y., Shibata, K., Hongo, A., *et al.* Ecabet sodium, a locally acting antiulcer drug, inhibits urease activity of *Helicobacter pylori*. *Eur. J. Pharmacol.* **345**(2), 193-198 (1998).
2. Hayashi, S., Sugiyama, T., Yachi, A., *et al.* Effect of ecabet sodium on *Helicobacter pylori* adhesion to gastric epithelial cells. *J. Gastroenterol.* **32**(5), 593-597 (1997).
3. Kinoshita, M., Iwasaki, H., Yasoshima, A., *et al.* Effects of ecabet sodium (TA-2711), a new antiulcer agent, on gastrointestinal mucosal prostanoid production and morphology in rats. *Biol. Pharm. Bull.* **16**(12), 1220-1225 (1993).
4. Okuyama, K., Saito, N., Kume, E., *et al.* Ecabet sodium prevents esophageal lesions induced by the reflux of gastric juice in rats. *Inflammopharmacology* **15**(2), 90-94 (2007).

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, 12/01/2022

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