

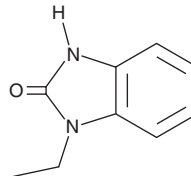
PRODUCT INFORMATION



1-EBIO

Item No. 14575

CAS Registry No.: 10045-45-1
Formal Name: 1-ethyl-1,3-dihydro-2H-benzimidazol-2-one
Synonym: 1-Ethylbenzimidazolinone
MF: C₉H₁₀N₂O
FW: 162.2
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 285 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

1-EBIO is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-EBIO in the solvent of choice, which should be purged with an inert gas. 1-EBIO is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1-EBIO in ethanol is approximately 30 mg/ml and approximately 100 mg/ml in DMSO and DMF.

1-EBIO is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1-EBIO should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 1-EBIO 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

1-EBIO is an activator of Ca²⁺-sensitive K⁺ channels (EC₅₀ = 490 μM in T84 human carcinoma cells).¹ This action is sensitive to the neurotoxin charybdotoxin (K_i = 3.5 nM). In mouse colonic epithelia, 1-EBIO also activates cAMP-sensitive K⁺ channels, a response this is inhibited by 293B.² In this type of epithelium, but not mouse nasal epithelia, 1-EBIO activates both types of channels, resulting in large Cl⁻ secretory currents.³ It also activates the human intermediate conductance Ca²⁺-activated K⁺ channel (EC₅₀ = 74 μM).⁴ 1-EBIO is currently used to study the role of these types of K⁺ channels in diverse physiological functions.^{5,6}

References

1. Devor, D.C., Singh, A.K., Frizzell, R.A., *et al.* Modulation of Cl⁻ secretion by benzimidazolones. I. Direct activation of a Ca(2+)-dependent K⁺ channel. *Am. J. Physiol.* **271**(5 Pt 1), L775-L784 (1996).
2. Cuthbert, A.W., Hickman, M.E., Thorn, P., *et al.* Activation of Ca²⁺- and cAMP-sensitive K⁺ channels in murine colonic epithelia by 1-ethyl-2-benzimidazolone. *Am. J. Physiol.* **277**(1), C111-C120 (1999).
3. MacVinish, L.J., Hickman, M.E., Mufti, D.A., *et al.* Importance of basolateral K⁺ conductance in maintaining Cl⁻ secretion in murine nasal and colonic epithelia. *J. Physiol.* **510**(1), 237-247 (1998).
4. Jensen, B.S., Strobaek, D., Christophersen, P., *et al.* Characterization of the cloned human intermediate-conductance Ca²⁺-activated K⁺ channel. *Am. J. Physiol.* **275**(3), C848-C856 (1998).
5. Gao, Z., Todorov, B., Barrett, C.F., *et al.* Cerebellar ataxia by enhanced Ca_v2.1 currents is alleviated by Ca²⁺-dependent K⁺-channel activators in Cacna1aS218L mutant mice. *J. Neurosci.* **32**(44), 15533-15546 (2012).
6. Maiz, J., Karakossian, M.H., Pakaprot, N., *et al.* Prolonging the postcomplex spike pause speeds eyeblink conditioning. *Proc. Natl. Acad. Sci. USA* **109**(41), 16726-16730 (2012).

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, 11/23/2022

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