

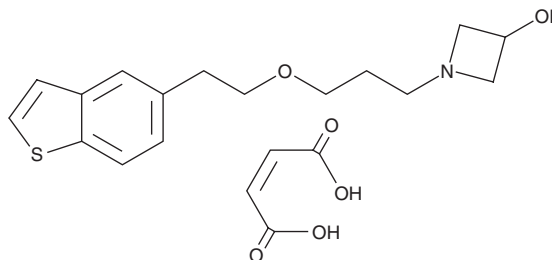
PRODUCT INFORMATION



Edonerpip (maleate)

Item No. 36026

CAS Registry No.: 519187-97-4
Formal Name: 1-[3-(2-benzo[b]thien-5-ylethoxy)propyl]-3-azetidinol, 2Z-butenedioate
Synonym: T-187MA
MF: $C_{16}H_{21}NO_2S \cdot C_4H_4O_4$
FW: 407.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 229 nm
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

Edonerpip (maleate) is supplied as a solid. A stock solution may be made by dissolving the edonerpip (maleate) in the solvent of choice, which should be purged with an inert gas. Edonerpip (maleate) is soluble in DMSO.

Description

Edonerpip is a neurotrophic agent.¹ It protects against sodium nitroprusside-induced mitochondrial dysfunction and apoptosis in primary rat cortical neurons when used at concentrations of 0.1 and 1 μM . Edonerpip (8.4 mg/kg per day, p.o.) prevents granule cell loss induced by amyloid- β (1-40) ($\text{A}\beta 40$) and improves performance in the distance movement, random reward place search, and place learning tasks in rats.² It also increases hippocampal neurogenesis and ameliorates $\text{A}\beta 40$ -induced spatial learning impairments in a rat model of Alzheimer's disease.³

References

1. Fukushima, T., Koide, M., Ago, Y., *et al.* T-817MA, a novel neurotrophic agent, improves sodium nitroprusside-induced mitochondrial dysfunction in cortical neurons. *Neurochem. Int.* **48**(2), 124-130 (2006).
2. Nguyen, P.T., Kimura, T., Ho, S.A., *et al.* Ameliorative effects of a neuroprotective agent, T-817MA, on place learning deficits induced by continuous infusion of amyloid-beta peptide (1-40) in rats. *Hippocampus* **17**(6), 443-455 (2007).
3. Kimura, T., Nguyen, P.T.H., Ho, S.A., *et al.* T-817MA, a neurotrophic agent, ameliorates the deficits in adult neurogenesis and spatial memory in rats infused i.c.v. with amyloid- β peptide. *Br. J. Pharmacol.* **157**(3), 451-463 (2009).

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, 12/19/2022

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