

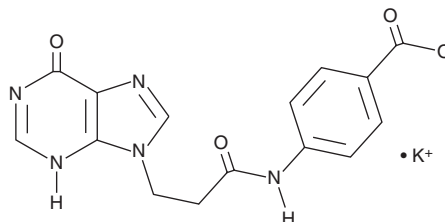
PRODUCT INFORMATION



Leteprinin (potassium salt)

Item No. 21748

CAS Registry No.: 192564-13-9
Formal Name: 4-[[3-(1,6-dihydro-6-oxo-9H-purin-9-yl)-1-oxopropyl]amino]-benzoic acid, monopotassium salt
Synonym: AIT 082
MF: C₁₅H₁₂N₅O₄ • K
FW: 365.4
Purity: ≥98%
UV/Vis.: λ_{max}: 265 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

Leteprinin (potassium salt) is supplied as a crystalline solid. Aqueous solutions of leteprinin (potassium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of leteprinin (potassium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Leteprinin is a nootropic and neuroprotective agent and a derivative of hypoxanthine (Item No. 22254).¹⁻³ Leteprinin (10 and 100 μM) enhances nerve growth factor induced neurite growth in PC12 cells and increases the production of synaptophysin when used at concentrations ranging from 5 to 100 ng/ml.^{1,3} It enhances working memory in a win-shift T-maze test in young mice and reduces time delay-induced working memory errors in old mice with mild or moderate, but not severe, age-induced memory deficits when administered at a dose of 30 mg/kg.⁴ In a rat model of spinal crush injury, leteprinin (60 mg/kg per day for 21 days) reduces impairments in locomotor function and reduces increases in the amount of necrotic tissue and the number of reactive astrocytes in the spinal cord.² It reduces hypoxia-induced decreases in neuronal density in the hippocampal CA1, CA3, and dentate gyrus regions and decreases the number of caspase-3 and TUNEL-positive cells in the hippocampus and prefrontal and parietal cortices in neonatal rats when administered at a dose of 60 mg/kg.⁵

References

1. Middlemiss, P.J., Glasky, A.J., Rathbone, M.P., *et al.* AIT-082, a unique purine derivative, enhances nerve growth factor mediated neurite outgrowth from PC12 cells. *Neurosci. Lett.* **199(2)**, 131-134 (1995).
2. Jiang, S., Khan, M.I., Middlemiss, P.J., *et al.* AIT-082 and methylprednisolone singly, but not in combination, enhance functional and histological improvement after acute spinal cord injury in rats. *Int. J. Immunopathol. Pharmacol.* **17(3)**, 353-366 (2004).
3. Lahiri, D.K., Ge, Y.-W., and Farlow, M.R. Effect of a memory-enhancing drug, AIT-082, on the level of synaptophysin. *Ann. N.Y. Acad. Sci.* **903(1)**, 387-393 (2000).
4. Glasky, A.J., Melchior, C.L., Pirzadeh, B., *et al.* Effect of AIT-082, a purine analog, on working memory in normal and aged mice. *Pharmacol. Biochem. Behav.* **47(2)**, 325-329 (1994).
5. Gencpinar, P., Tüzün, F., Ozbal, S., *et al.* Effects of neotrofin on neonatal hypoxic ischemic brain injury. *Neurosci. Lett.* **505(2)**, 205-210 (2011).

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, 11/07/2022

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