

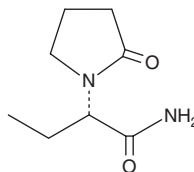
PRODUCT INFORMATION



Levetiracetam

Item No. 9001820

CAS Registry No.: 102767-28-2
Formal Name: (αS)-ethyl-2-oxo-1-pyrrolidineacetamide
Synonym: UCB-L 059
MF: C₈H₁₄N₂O₂
FW: 170.2
Purity: ≥98%
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

Levetiracetam is supplied as a crystalline solid. A stock solution may be made by dissolving the levetiracetam in the solvent of choice. Levetiracetam is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of levetiracetam in these solvents is approximately 16.7, 20, and 20 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 levetiracetam can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of levetiracetam in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Levetiracetam is an antiepileptic agent that binds to the synaptic vesicle protein SV2A (IC₅₀ = 1.99 μM).¹ *In vivo*, levetiracetam suppresses seizures in fully kindled rats.²

References

1. Lynch, B.A., Lamberg, N., Nocka, K., et al. *Proc. Natl. Acad. Sci. USA* **101**(26), 9861-9866 (2004).
2. Löscher, W. and Brandt, C. *Pharmacol. Rev.* **62**(4), 668-700 (2010).

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, 11/21/2022

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