

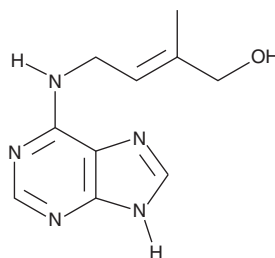
PRODUCT INFORMATION



trans-Zeatin

Item No. 13226

CAS Registry No.: 1637-39-4
Formal Name: 2E-methyl-4-(9H-purin-6-ylamino)-2-buten-1-ol
MF: C₁₀H₁₃N₅O
FW: 219.2
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 269 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

trans-Zeatin is supplied as a crystalline solid. A stock solution may be made by dissolving the *trans*-zeatin in the solvent of choice, which should be purged with an inert gas. *trans*-Zeatin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of *trans*-zeatin in these solvents is approximately 3 and 0.5 mg/ml, respectively.

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

trans-Zeatin is a cytokinin plant growth regulator with antioxidant and neuroprotective activities.¹⁻⁴ It binds to the cytokinin receptor *Arabidopsis* histidine kinase 3 (AHK3) with a K_D value of 1.3 nM.⁵ *trans*-Zeatin increases chlorophyll levels in etiolated *Cucumis sativus* cotyledons in a concentration-dependent manner.¹ It increases callus growth and shoot formation in *N. tabacum* calluses when used at concentrations of 5 and 50 μM.² *trans*-Zeatin (25-100 μM) reduces production of reactive oxygen species (ROS) induced by amyloid β (25-35) (Aβ₂₅₋₃₅) in PC12 cells.⁴ It reduces scopolamine-induced spontaneous alternations in the Y-maze, indicating enhanced spatial memory, in mice when administered at doses of 1.5, 3, and 4.5 mg/kg per day.

References

1. Fletcher, R.A. and McCullagh, D. Cytokinin-induced chlorophyll formation in cucumber cotyledons. *Planta* **101**(1), 88-90 (1971).
2. Yamada, Y., Sekiya, J., and Koshimizu, K. Cytokinin-induced shoot formation. *Phytochemistry* **11**(3), 1019-1021 (1972).
3. Romanov, G.A., Lomin, S.N., and Schmölling, T. Biochemical characteristics and ligand-binding properties of *Arabidopsis* cytokinin receptor AHK3 compared to CRE1/AHK4 as revealed by a direct binding assay. *J. Exp. Bot.* **57**(15), 4051-4058 (2006).
4. Choi, S.J., Jeong, C.-H., Choi, S.-G., et al. Zeatin prevents amyloid β-induced neurotoxicity and scopolamine-induced cognitive deficits. *J. Med. Food* **12**(2), 271-277 (2009).
5. Romanov, G.A., Lomin, S.N., and Schmölling, T. Biochemical characteristics and ligand-binding properties of *Arabidopsis* cytokinin receptor AHK3 compared to CRE1/AHK4 as revealed by a direct binding assay. *J. Exp. Bot.* **57**(15), 4051-4058 (2006).

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, 12/13/2022

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