

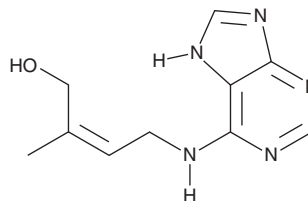
# PRODUCT INFORMATION



## cis-Zeatin

Item No. 35549

CAS Registry No.: 32771-64-5  
Formal Name: 2Z-methyl-4-(9H-purin-6-ylamino)-2-buten-1-ol  
MF:  $C_{10}H_{13}N_5O$   
FW: 219.2  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

*cis*-Zeatin is supplied as a solid. A stock solution may be made by dissolving the *cis*-zeatin in the solvent of choice, which should be purged with an inert gas. *cis*-Zeatin is soluble in methanol and DMSO.

### Description

*cis*-Zeatin is a cytokinin plant growth regulator and an isomer of *trans*-zeatin (Item No. 13226).<sup>1</sup> It induces callus growth in isolated *N. tabacum* callus tissue when used at concentrations ranging from 0.1 to 1  $\mu\text{M}$ .

### Reference

1. Schmitz, R.Y. and Skoog, F. Cytokinins: Synthesis and biological activity of geometric and position isomers of zeatin. *Plant Physiol.* **50**(6), 702-705 (1972).

#### 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, 12/12/2022

#### 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