

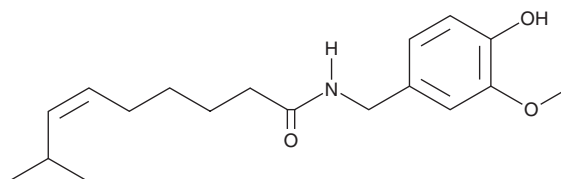
# PRODUCT INFORMATION



## (Z)-Capsaicin

Item No. 39929

**CAS Registry No.:** 25775-90-0  
**Formal Name:** N-[(4-hydroxy-3-methoxyphenyl)methyl]-8-methyl-6-nonenamide  
**Synonyms:** *cis*-Capsaicin, *cis*-8-methyl-N-Vanillyl-6-nonenamide, Zucapsaicin  
**MF:** C<sub>18</sub>H<sub>27</sub>NO<sub>3</sub>  
**FW:** 305.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

(Z)-Capsaicin is supplied as a solid. A stock solution may be made by dissolving the (Z)-capsaicin in the solvent of choice, which should be purged with an inert gas. (Z)-Capsaicin is soluble in methanol and DMSO.

### Description

(Z)-Capsaicin is an alkaloid that has been found in *Capsicum*, has diverse biological activities, and is the *cis*-isomer of capsaicin (Item Nos. 92350 | 10010743).<sup>1-6</sup> It is an agonist of transient receptor potential vanilloid 1 (TRPV1), inducing calcium uptake in CHO cells expressing human TRPV1 (EC<sub>50</sub> = 28.2 nM).<sup>2</sup> (Z)-Capsaicin (10 μM) shortens the action potential duration in isolated canine Purkinje fibers.<sup>3</sup> It increases the latency to paw withdrawal in the hot plate test in rats when administered at a dose of 0.68 mmol/animal.<sup>4</sup> Topical application of (Z)-capsaicin (25 μl of a 1% solution) reduces croton oil-induced ear edema in mice. It induces emesis in musk shrews (*S. murinus*).<sup>5</sup> (Z)-Capsaicin decreases disease severity in a guinea pig model of genital herpes induced by herpes simplex virus 2 (HSV-2).<sup>6</sup> Formulations containing (Z)-capsaicin have been used in the treatment of osteoarthritis and neuropathic pain.

### References

1. Espichán, F., Rojas, R., Quispe, F., *et al.* Metabolomic characterization of 5 native Peruvian chili peppers (*Capsicum* spp.) as a tool for species discrimination. *Food Chem.* **386**, 132704 (2022).
2. Ann, J., Kim, H.S., Thorat, S.A., *et al.* Discovery of nonpungent transient receptor potential vanilloid 1 (TRPV1) agonist as strong topical analgesic. *J. Med. Chem.* **63**(1), 418-424 (2020).
3. Arnar, D.O., Cai, J.J., Lee, H.-C., *et al.* Electrophysiologic effects of civamide (zucapsaicin) on canine cardiac tissue in vivo and in vitro. *J. Cardiovasc. Pharmacol.* **32**(6), 875-883 (1998).
4. Janusz, J.M., Buckwalter, B.L., Young, P.A., *et al.* Vanilloids. 1. Analogs of capsaicin with antinociceptive and antiinflammatory activity. *J. Med. Chem.* **36**, 2595-2604 (1993).
5. Rudd, J.A. and Wai, M.K. Genital grooming and emesis induced by vanilloids in *Suncus murinus*, the house musk shrew. *Eur. J. Pharmacol.* **422**(1-3), 185-195 (2001).
6. Bourne, N., Bernstein, D.I., and Stanberry, L.R. Civamide (*cis*-capsaicin) for treatment of primary or recurrent experimental genital herpes. *Antimicrob. Agents Chemother.* **43**(11), 2685-2688 (1999).

#### 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, 02/19/2024

#### 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