

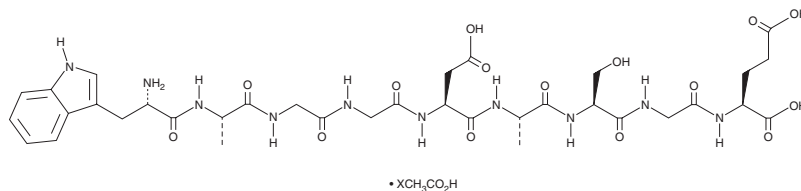
# PRODUCT INFORMATION



## Emideltide (acetate)

Item No. 34426

<b>Synonyms:</b>	Delta Sleep-inducing Peptide, DSIP
<b>MF:</b>	$C_{35}H_{48}N_{10}O_{15} \cdot XC_2H_4O_2$
<b>FW:</b>	848.8
<b>Purity:</b>	≥98%
<b>UV/Vis.:</b>	$\lambda_{max}$ : 221 nm
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years



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

### Laboratory Procedures

Emideltide (acetate) is supplied as a solid. A stock solution may be made by dissolving the emideltide (acetate) in the solvent of choice, which should be purged with an inert gas. Emideltide (acetate) is slightly soluble in DMSO.

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 emideltide (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of emideltide (acetate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Emideltide, also known as delta sleep-inducing peptide (DSIP), is a synthetic nonapeptide sedative.<sup>1</sup> *In vivo*, emideltide (30 nmol/kg, i.v.) increases EEG delta activity and induces orthodox slow-wave sleep in free-moving rabbits. It also increases non-rapid eye movement (NREM) and total sleep in free-moving rats.<sup>2</sup>

### References

1. Monnier, M., Dudler, L., Gächter, R., *et al.* Delta sleep-inducing peptide (DSIP): EEG and motor activity in rabbits following intravenous administration. *Neurosci. Lett.* **6(1)**, 9-13 (1977).
2. Kafi, S., Monnier, M., and Gallard, J.M. The delta-sleep inducing peptide (DSIP) increases duration of sleep in rats. *Neurosci. Lett.* **13(2)**, 169-172 (1979).

#### 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, 11/22/2022

#### 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