

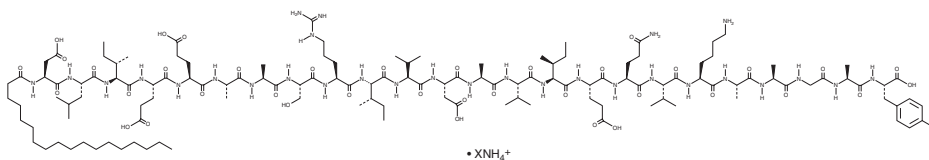
PRODUCT INFORMATION



st-Ht31 (ammonium salt)

Item No. 36745

Synonym: sHt31
Peptide Sequence: stearoyl-DLIEEAASRIV
DAVIEQVKAAGAY-OH
MF: C₁₂₉H₂₁₇N₂₉O₃₉ • XNH₄
FW: 2,798.3
Purity: ≥98%
Supplied as: A 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

st-Ht31 (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the st-Ht31 (ammonium salt) in the solvent of choice, which should be purged with an inert gas. st-Ht31 (ammonium salt) is soluble in 30% acetonitrile/water.

Description

st-Ht31 is a cell-permeable steared peptide derived from the amphipathic helix domain from A-kinase anchoring proteins (AKAPs), which binds protein kinase A (PKA).^{1,2} It decreases PKA membrane anchoring but not activity in a reporter assay using BHK-21 fibroblasts when used at a concentration of 50 μM.² st-Ht31 (1 and 3 μM) reduces capacitance and hyperactivation, as well as prevents *in vitro* fertilization, in mouse sperm.¹ It increases cholesterol export from BHK-21 cells and RAW 264.7 macrophages expressing ATP-binding cassette transporter 1 (ABCA1) when used at a concentration of 50 μM.²

References

1. Stival, C., Ritagliati, C., Xu, X., *et al.* Disruption of protein kinase A localization induces acrosomal exocytosis in capacitated mouse sperm. *J. Biol. Chem.* **293**(23), 9435-9447 (2018).
2. Ma, L., Dong, F., Denis, M., *et al.* Ht31, a protein kinase A anchoring inhibitor, induces robust cholesterol efflux and reverses macrophage foam cell formation through ATP-binding cassette transporter A1. *J. Biol. Chem.* **286**(5), 3370-3378 (2011).

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, 10/11/2023

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