

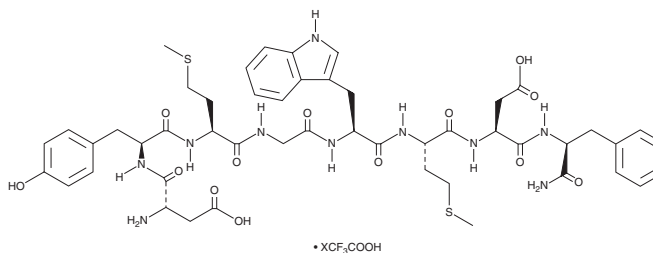
PRODUCT INFORMATION



CCK Octapeptide (non-sulfated) (trifluoroacetate salt)

Item No. 24404

Formal Name: 2-desulfo-cholecystokinin-8 (swine), trifluoroacetate salt
Synonyms: Cholecystokinin Octapeptide (26-33, Pancreozymin (C-terminal) Octapeptide, SQ 19,265)
MF: C₄₉H₆₂N₁₀O₁₃S₂ • XCF₃COOH
FW: 1,063.2
Purity: ≥95%
Supplied as: A lyophilized powder
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Cholecystokinin (CCK) octapeptide (non-sulfated) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the CCK octapeptide (non-sulfated) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. CCK octapeptide (non-sulfated) (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

CCK octapeptide is a peptide hormone found in the intestine and brain that stimulates digestion, mediates satiety, and is involved in anxiety.^{1,2} It is a cleavage product of the full CCK preprohormone. CCK octapeptide (non-sulfated) has 1/300th of the activity of sulfated CCK octapeptide.³

References

1. Raybould, H.E. and Lloud, K.C.K. Integration of postprandial function in the proximal gastrointestinal tract. Role of CCK and sensory pathways. *Ann. N.Y. Acad. Sci.* **713**, 143-156 (1994).
2. Lydiard, R.B. Neuropeptides and anxiety: Focus on cholecystokinin. *Clin. Chem.* **40(2)**, 315-318 (1994).
3. Ondetti, M.A., Plušćec, J., Sabo, E.F., *et al.* Synthesis of cholecystokinin-pancreozymin. I. The C-terminal dodecapeptide. *J. Am. Chem. Soc.* **92(1)**, 195-199 (1970).

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, 12/20/2022

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