

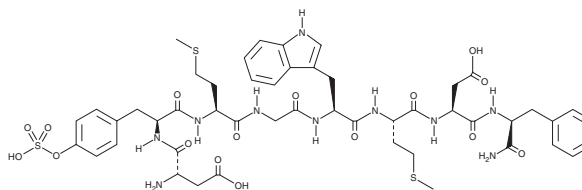
PRODUCT INFORMATION



CCK Octapeptide (sulfated)

Item No. 23371

CAS Registry No.: 25126-32-3
Formal Name: 5-L-methionine-3-10-caerulein
Synonyms: CCK-8S, Cholecystokinin (26-33), SQ 19,844, Sincalide
MF: C₄₉H₆₂N₁₀O₁₆S₃
FW: 1,143.3
Purity: ≥98%
Supplied as: A crystalline 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

CCK octapeptide (sulfated) is supplied as a crystalline solid. Aqueous solutions of CCK octapeptide (sulfated) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CCK octapeptide (sulfated) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cholecystokinin (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 cleaved product of the full CCK prohormone that is more potent than the full length enzyme in stimulating contraction of guinea pig gallbladder but with a shorter duration.^{2,3} CCK octapeptide (sulfated) is 300-fold more active than the non-sulfated CCK octapeptide (Item No. 24404).²

References

1. Raybould, H.E. and Lloyd, K.C. Integration of postprandial function in the proximal gastrointestinal tract. Role of CCK and sensory pathways. *Ann. N.Y. Acad. Sci.* **713(1)**, 143-156 (1994).
2. Ondetti, M.A., Plušćec, J., and Sabo, E.F. Synthesis of cholecystokinin-pancreozymin. I. The C-terminal dodecapeptide. *J. Am. Chem. Soc.* **92(1)**, 195-199 (1970).
3. Rubin, B., Engel, S.L., Drungis, A.M., *et al.* Cholecystokinin-like activities in guinea pigs and in dogs of the C-terminal octapeptide (SQ 19,844) of cholecystokinin. *J. Pharm. Sci.* **58(8)**, 955-959 (1969).

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, 12/13/2022

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