

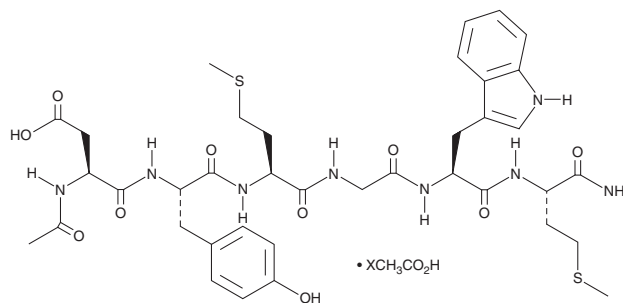
# PRODUCT INFORMATION



## CCK (26-31) (non-sulfated) (acetate)

Item No. 24951

**Formal Name:** N-acetyl-L- $\alpha$ -aspartyl-L-tyrosyl-L-methionylglycyl-L-tryptophyl-L-methioninamide, acetate  
**Synonym:** N-Acetyl Cholecystokinin (26-31), (non-sulfated)  
**MF:** C<sub>38</sub>H<sub>50</sub>N<sub>8</sub>O<sub>10</sub>S<sub>2</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>  
**FW:** 843.0  
**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

CCK (26-31) (non-sulfated) (acetate) is supplied as a lyophilized powder. A stock solution may be made by dissolving the CCK (26-31) (non-sulfated) (acetate) in the solvent of choice, which should be purged with an inert gas. CCK (26-31) (non-sulfated) (acetate) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

### Description

CCK (26-31) is an N-terminal fragment of CCK (Item Nos. 23371 | 24404), a peptide hormone found in the intestine and brain that stimulates digestion, mediates satiety, and is involved in anxiety.<sup>1,2</sup>

### 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., 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, 03/06/2026

#### 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