

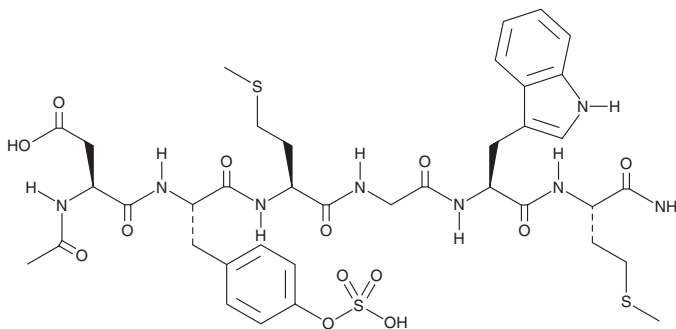
# PRODUCT INFORMATION



## CCK (26-31) (sulfated)

Item No. 24950

**CAS Registry No.:** 89911-65-9  
**Formal Name:** N-acetyl-L- $\alpha$ -aspartyl-O-sulfo-L-tyrosyl-L-methionylglycyl-L-tryptophyl-L-methioninamide  
**Synonym:** N-Acetyl Cholecystokinin (26-31), (sulfated)  
**MF:** C<sub>38</sub>H<sub>50</sub>N<sub>8</sub>O<sub>13</sub>S<sub>3</sub>  
**FW:** 923.0  
**Purity:**  $\geq 95\%$   
**Supplied as:** A lyophilized powder  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

CCK (26-31) (sulfated) is supplied as a lyophilized powder. A stock solution may be made by dissolving the CCK (26-31) (sulfated) in water. The solubility of CCK (26-31) (sulfated) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### 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> The sulfated form of CCK (26-31) inhibits binding of [<sup>125</sup>I]CCK-33 to guinea pig cortical membranes by 21% when used at a concentration of 0.1 mM.<sup>3</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., Pluščec, J., Sabo, E.F., *et al.* Synthesis of cholecystokinin-pancreozymin. I. The C-terminal dodecapeptide. *J. Am. Chem. Soc.* **92**(1), 195-199 (1970).
3. Knight, M., Tamminga, C.A., Steardo, L., *et al.* Cholecystokinin-octapeptide fragments: Binding to brain cholecystokinin receptors. *Eur. J. Pharmacol.* **105**(1-2), 49-55 (1984).

#### 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, 11/18/2022

#### 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