

# PRODUCT INFORMATION



## Biocytin-β-Endorphin (human) (trifluoroacetate salt)

Item No. 25011

**Formal Name:** N-[N<sup>6</sup>-[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]-L-lysine]-27-L-tyrosine-31-L-glutamic acid-β-endorphin, trifluoroacetate salt

**MF:** C<sub>174</sub>H<sub>277</sub>N<sub>43</sub>O<sub>49</sub>S<sub>2</sub> • XCF<sub>3</sub>COOH

**FW:** 3,819.5

**Purity:** ≥95%

**Supplied as:** A lyophilized powder

**Storage:** -20°C

**Stability:** ≥4 years

Biotin—Lys—Tyr—Gly—Gly—Phe—Met—Thr—Ser—Glu—Lys—Ser—  
Gln—Thr—Pro—Leu—Val—Thr—Leu—Phe—Lys—Asn—  
Ala—Ile—Ile—Lys—Asn—Ala—Tyr—Lys—Lys—Gly—Glu—OH

• XCF<sub>3</sub>COOH

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

### Laboratory Procedures

Biocytin-β-endorphin (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the biocytin-β-endorphin (human) (trifluoroacetate salt) in water. The solubility of biocytin-β-endorphin (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Biocytin-β-endorphin is a form of β-endorphin (Item Nos. 24955 | 24153) that is conjugated to biocytin. It has been used to examine β-endorphin localization and binding site dynamics in live *R. esculenta* pars distalis pituitary cells.<sup>1</sup>

### Reference

1. Sabbieti, M.G., Marchetti, L., Menghi, G., *et al.* Occurrence of β-endorphin binding sites in the pituitary of the frog *Rana esculenta*: Effect of β-endorphin on luteinizing hormone secretion. *Gen. Comp. Endocrinol.* **132**(3), 391-398 (2003).

#### 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, 11/21/2022

#### 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