

PRODUCT INFORMATION



Tyr-β-CGRP (human) (trifluoroacetate salt)

Item No. 24766

Synonym: Tyr-β-Calcitonin Gene-Related Peptide (human), [Tyr⁰] Calcitonin Gene-Related Peptide-2 (human), Tyr-CGRP-2 (human)

MF: C₁₇₁H₂₇₆N₅₂O₅₀S₃ • XCF₃COOH

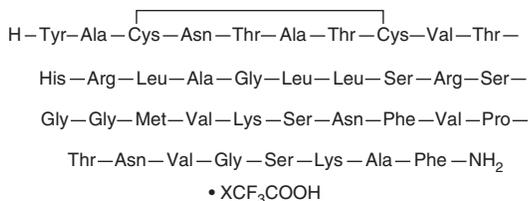
FW: 3,956.5

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

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

Description

Tyr-β-CGRP is an N-terminal extended tyrosinated analog of β-calcitonin gene-related peptide (β-CGRP; Item No. 24725).

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