

PRODUCT INFORMATION



Katacalcin (trifluoroacetate salt)

Item No. 39285

Formal Name: L- α -aspartyl-L-methionyl-L-seryl-L-seryl-L- α -aspartyl-L-leucyl-L- α -glutamyl-L-arginyl-L- α -aspartyl-L-histidyl-L-arginyl-L-prolyl-L-histidyl-L-valyl-L-seryl-L-methionyl-L-prolyl-L-glutamyl-L-asparaginyll-L-alanyl-L-asparagine, trifluoroacetate salt
H—Asp—Met—Ser—Ser—Asp—Leu—Glu—Arg—Asp—His—
Arg—Pro—His—Val—Ser—Met—Pro—Gln—Asn—Ala—
Asn—OH
• XCF₃COOH

Synonym: PDN-21

Peptide Sequence: DMSSDLERDHRPHVSMQPQAN-OH

MF: C₉₇H₁₅₄N₃₄O₃₆S₂ • XCF₃COOH

FW: 2,436.6

Purity: ≥95%

Supplied as: A 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

Katacalcin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the katacalcin (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Katacalcin (trifluoroacetate salt) is soluble in methanol. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Katacalcin is an endogenous peptide (121-141) encoded by the CALCA gene and is expressed in, and secreted from, thyroid parafollicular cells in humans.¹ Unlike calcitonin (Item No. 24409), a calcium regulating peptide hormone, katacalcin does not affect serum calcium levels. Plasma levels of katacalcin are increased in patients with medullary thyroid carcinoma.²

References

1. Fischer, J.A. and Born, W. Novel peptides from the calcitonin gene: Expression, receptors and biological function. *Peptides* **6(Suppl 3)**, 265-271 (1985).
2. Takami, H., Shikata, J.-I., Horie, H., et al. PDN-21: Possible tumor marker for medullary thyroid carcinoma. *J. Surg. Oncol.* **44(4)**, 205-207 (1990).

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, 09/26/2023

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