

PRODUCT INFORMATION



Parathyroid Hormone (1-13) (trifluoroacetate salt)

Item No. 24984

Formal Name:	L-seryl-L-valyl-L-seryl-L- α -glutamyl-L-isoleucyl-L-glutamyl-L-leucyl-L-methionyl-L-histidyl-L-asparaginyl-L-leucylglycyl-L-lysine, trifluoroacetate salt	H — Ser — Val — Ser — Glu — Ile — Gln — Leu — Met — His — Asn —
Synonym:	PTH (1-13)	Leu — Gly — Lys — OH
MF:	$C_{62}H_{106}N_{18}O_{20}S \cdot XCF_3COOH$	$\bullet XCF_3COOH$
FW:	1,455.7	
Purity:	$\geq 95\%$	
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

Parathyroid hormone (PTH) (1-13) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the PTH (1-13) (trifluoroacetate salt) in water. The solubility of PTH (1-13) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PTH (1-13) is an N-terminal fragment of PTH (1-34) (Item No. 24985).¹ *In vivo*, PTH (1-13) (0.003-1 μ mol/kg) increases mean arterial pressure (MAP) in anesthetized rats.

Reference

1. Nakamura, R., Sokabe, H., Kimura, T., *et al.* Action of fragments of human parathyroid hormone on blood pressure in rats. *Endocrinol. Jpn.* **28**(4), 547-549 (1981).

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, 12/19/2022

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