

PRODUCT INFORMATION

Orexin A (17-33) (human, mouse, rat, bovine) (trifluoroacetate salt) Item No. 36765

Formal Name:	L-tyrosyl-L- α -glutamyl-L-leucyl-L-leucyl-L-histidylglycyl-L-alanylglycyl-L-asparaginy-L-histidyl-L-alanyl-L-alanylglycyl-L-isoleucyl-L-leucyl-L-threonyl-L-leucinamide, trifluoroacetate salt	H-Tyr-Glu-Leu-Leu-His-Gly-Ala-Gly-Asn-His-Ala-Ala-Gly-Ile-Leu-Thr-Leu-NH ₂
Synonym:	OXA (17-33)	
MF:	C ₇₉ H ₁₂₅ N ₂₃ O ₂₂ • XCF ₃ COOH	
FW:	1,749.0	• XCF ₃ COOH
Purity:	≥98%	
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

Orexin A (17-33) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the orexin A (17-33) (human, mouse, rat, bovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Orexin A (17-33) (human, mouse, rat, bovine) (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of orexin A (17-33) (human, mouse, rat, bovine) (trifluoroacetate salt) in these solvents is approximately 5 mg/ml and 1 mg/ml, respectively.

Description

Orexin A (17-33) is a truncated form of orexin A and peptide orexin receptor 1 (OX1R) agonist.¹ It selectively induces calcium mobilization in CHO cells overexpressing human OX1R over CHO cells overexpressing human OX2R (EC₅₀s = 8.29 and 187 nM, respectively).

Reference

- German, N.A., Decker, A.M., Gilmour, B.P., *et al.* Truncated orexin peptides: Structure-activity relationship studies. *ACS Med. Chem. Lett.* **4**(12), 1224-1227 (2013).

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, 12/21/2022

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