

PRODUCT INFORMATION

FQWQRNIRKVR-NH₂ (trifluoroacetate salt)

Item No. 42400

Formal Name:	L-phenylalanyl-L-glutamyl-L-tryptophyl-L-glutamyl-L-arginyl-L-asparagyl-L-isoleucyl-L-arginyl-L-lysyl-L-valyl-L-argininamide, trifluoroacetate salt	H—Phe—Gln—Trp—Gln—Arg—Asn—Ile—Arg—Lys—Val—
Synonym:	LF11	Arg—NH ₂
Peptide Sequence:	FQWQRNIRKVR-NH ₂	
MF:	C ₆₉ H ₁₁₂ N ₂₆ O ₁₄ • XCF ₃ COOH	• XCF ₃ COOH
FW:	1,529.8	
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

FQWQRNIRKVR-NH₂ (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the FQWQRNIRKVR-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. FQWQRNIRKVR-NH₂ (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Description

FQWQRNIRKVR-NH₂ is a peptide fragment of lactoferricin with antibacterial activity.¹ It is active against *E. coli* (MIC = 256 µg/ml).

References

1. Zweytick, D., Pabst, G., Abuja, P.M., *et al.* Influence of N-acylation of a peptide derived from human lactoferricin on membrane selectivity. *Biochim. Biophys. Acta* **1758(9)**, 1426-1435 (2006).

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, 01/24/2025

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