

PRODUCT INFORMATION



(DYKDDDDK)₃ (trifluoroacetate salt)

Item No. 44482

Peptide Sequence: DYKDDDDKDYKDDDDKDYKDDDDK-OH

MF: C₁₂₃H₁₇₆N₃₀O₅₈ • XCF₃COOH

FW: 3,002.9

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H-Asp-Tyr-Lys-Asp-Asp-Asp-Asp-Lys-Asp-Tyr-

Lys-Asp-Asp-Asp-Asp-Lys-Asp-Tyr-Lys-Asp-

Asp-Asp-Asp-Lys-OH

• XCF₃COOH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(DYKDDDDK)₃ (trifluoroacetate salt) is supplied as a solid. Aqueous solutions of (DYKDDDDK)₃ (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. (DYKDDDDK)₃ (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

(DYKDDDDK)₃ is a peptide containing three tandem repeats of the epitope tag DYKDDDDK that has been used to elute DYKDDDDK-tagged proteins during affinity purification.¹

Reference

1. Han, Z. and Svejstrup, J.Q. Protocol to detect spontaneous termination of yeast RNAPII transcription *in vitro*. *STAR Protoc.* 5(4), 103369 (2024).

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, 12/03/2025

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