

PRODUCT INFORMATION



POLA2 (22-30) (human) (trifluoroacetate salt)

Item No. 44402

CAS Registry No.: 1000281-91-3
Formal Name: L-alanyl-L-leucyl-L-isoleucyl-L- α -glutamyl-L-lysyl-L-leucyl-L-valyl-L- α -glutamyl-L-leucine
Synonyms: Ala-Leu-Ile-Glu-Lys-Leu-Val-Glu-Leu-OH,
DNA Polymerase α Subunit B (22-30), H-Ala-Leu-Ile-Glu-Lys-Leu-Val-Glu-Leu-OH
p70 (22-30)
Peptide Sequence: ALIEKLVEL-OH • XCF₃COOH
MF: C₄₈H₈₆N₁₀O₁₄ • XCF₃COOH
FW: 1,027.3
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years

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

Laboratory Procedures

POLA2 (22-30) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the POLA2 (22-30) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. POLA2 (22-30) (human) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO.

Description

DNA polymerase α subunit B (POLA2) (22-30) is a peptide fragment of POLA2, which is a protein member of DNA polymerase α and involved in initiating DNA replication and DNA double-strand break (DSB) repair.

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, 10/21/2025

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