

PRODUCT INFORMATION



DNAJ-PKAc Fusion Peptide (24-mer) (trifluoroacetate salt)

Item No. 45191

Formal Name: (S)-N1-((S)-1-(((3S,4S,7S,12S)-12-(2-((5S,8S)-1,8-diamino-11-((diaminomethylene)amino)-6,7-dioxoundecan-5-yl)hydrazineyl)-15-((diaminomethylene)amino)-7-(hydroxyamino)-3-methyl-5,6,10,11-tetraoxopentadecan-4-yl)amino)-1-oxo-3-phenylpropan-2-yl)-N4-((6S,9S,17S,22S,25S,28S,33S,36S,39S,42S,45S,48S,51S,56S,59S,63S)-1,1-diamino-28,45,51-tris(4-aminobutyl)-36-benzyl-17,22,33,56,59,63-hexakis(hydroxyamino)-9-(4-hydroxybenzyl)-39-isobutyl-25-isopropyl-42,48-dimethyl-7,8,11,14,18,19,23,24,27,30,34,35,38,41,44,-47,50,53,57,58,61,62-docosaono-64-phenyl-2,10,13,26,29,37,40,43,46,49,52-undecaazatetrahexacont-1-en-6-yl)-2-(hydroxyamino)succinamide, trifluoroacetate

H—Arg—Lys—Arg—Glu—Ile—Phe—Asp—Arg—Tyr—Gly—
Glu—Glu—Val—Lys—Glu—Phe—Leu—Ala—Lys—Ala—
Lys—Glu—Asp—Phe—OH
• XCF₃COOH

Synonyms: DNAJB1-PRKACA Fusion,
Hepatocellular Carcinoma 24-mer Fusion

Peptide Sequence: RKREIFDRYGEEVKEFLAKAKEDF-OH

MF: C₁₃₆H₂₁₁N₃₇O₄₀ • XCF₃COOH

FW: 3,004.4

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

DNAJ-PKAc fusion peptide (24-mer) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the DNAJ-PKAc fusion peptide (24-mer) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. DNAJ-PKAc fusion peptide (24-mer) (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of DNAJ-PKAc fusion peptide (24-mer) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. DNAJ-PKAc fusion peptide (24-mer) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

DNAJ-PKAc Fusion Peptide (24-mer) is a synthetic peptide spanning the junction of the DNAJ-PKAc fusion protein found in fibrolamellar carcinoma (FLC) tumors.¹ It contains amino acids 59-70 of DNAJ heat shock protein family member B1 (DNAJB1) and 16-27 of PKA α , which are located in the J domain and catalytic domain, respectively.

Reference

1. O'Neill, A.F., Zeme, E.L., Abou-Alfa, G.K., et al. *Hepatology* (2026).

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, 04/10/2026

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