

# PRODUCT INFORMATION



## KLQEFLQTL (trifluoroacetate salt)

Item No. 44401

CAS Registry No.: 474283-38-0

Formal Name: L-lysyl-L-leucyl-L-glutamyl-L- $\alpha$ -glutamyl-L-phenylalanyl-L-leucyl-L-glutamyl-L-threonyl-L-leucine, trifluoroacetate salt

Synonym: H-Lys-Leu-Gln-Glu-Phe-Leu-Gln-Thr-Leu-OH      H-Lys-Leu-Gln-Glu-Phe-Leu-Gln-Thr-Leu-OH  
• XCF<sub>3</sub>COOH

Peptide Sequence: H-KLQEFLQTL-OH

MF: C<sub>52</sub>H<sub>86</sub>N<sub>12</sub>O<sub>15</sub> • XCF<sub>3</sub>COOH

FW: 1,119.3

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

KLQEFLQTL (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the KLQEFLQTL (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. KLQEFLQTL (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO. KLQEFLQTL (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS pH 7.2.

### Description

KLQEFLQTL is a peptide corresponding to amino acids 16-24 of Fanconi anemia group I protein (FACI).

#### 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, 04/27/2026

#### 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