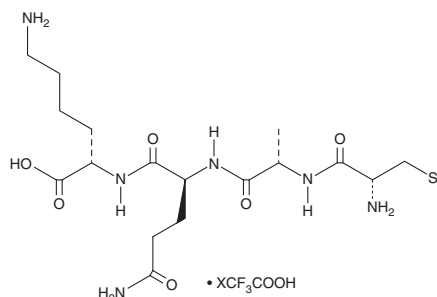


PRODUCT INFORMATION

CAQK Targeting/Delivery Peptide (trifluoroacetate salt)

Item No. 45529

Formal Name: L-cysteinyl-L-alanyl-L-glutaminyl-L-lysine
Synonyms: CAQK Peptide, H-Cys-Ala-Gln-Lys-OH
Peptide Sequence: H-CAQK-OH
MF: $C_{17}H_{32}N_6O_6S \cdot XCF_3COOH$
FW: 448.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

CAQK targeting/delivery peptide (CAQK peptide) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the CAQK peptide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. CAQK peptide (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in ethanol and 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 CAQK peptide (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. CAQK peptide (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

CAQK peptide is a four-amino acid synthetic peptide that interacts with versican and tenascin-R, proteoglycans that are upregulated in the brain extracellular matrix (ECM) in CNS injuries.¹ Intravenous administration of CAQK peptide leads to its accumulation in injured brain tissue over uninjured brain tissue in mice. CAQK peptide-conjugated nanoparticles have been used to selectively deliver siRNA to injured brain tissue and to image brain injuries in mice.

Reference

1. Mann, A.P., Scodeller, P., Hussain, S., et al. A peptide for targeted, systemic delivery of imaging and therapeutic compounds into acute brain injuries. *Nat. Commun.* **7**, 11980 (2016).

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, 06/03/2026

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