

PRODUCT INFORMATION

Caveolin-1 (82-101) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt)

Item No. 37003

Formal Name:	L- α -aspartylglycyl-L-isoleucyl-L-tryptophyl-L-lysyl-L-alanyl-L-seryl-L-phenylalanyl-L-threonyl-L-threonyl-L-phenylalanyl-L-threonyl-L-valyl-L-threonyl-L-lysyl-L-tyrosyl-L-tryptophyl-L-phenylalanyl-L-tyrosyl-L-arginine, trifluoroacetate salt	H—Asp—Gly—Ile—Trp—Lys—Ala—Ser—Phe—Thr—Thr—Phe—Thr—Val—Thr—Lys—Tyr—Trp—Phe—Tyr—Arg—NH ₂
Synonyms:	Cav-1, Caveolin 1 Scaffolding Domain Peptide, CSD, CSP, DGIWKASFTTFTVTKYWFYR-NH ₂	• XCF ₃ COOH
MF:	C ₁₂₄ H ₁₇₀ N ₂₈ O ₂₉ • XCF ₃ COOH	
FW:	2,516.9	
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

Caveolin-1 (82-101) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the caveolin-1 (82-101) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Caveolin-1 (82-101) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 3 mg/ml.

Description

Caveolin-1 (82-101) amide is a 20-amino acid synthetic peptide fragment corresponding to residues 82-101 of caveolin-1, which is involved in endocytosis and interacts with a variety of receptors and other proteins.¹ Caveolin-1 (82-101) (0.8 μ mol/kg) reduces aging-induced increases in microvascular leakage and fibrosis, as well as the ratio of phosphorylated-to-unphosphorylated PDGFR and protein-rich tyrosine 2 (Pyk2), in aged mouse heart, kidney, and brain. It also increases left ventricular ejection fraction and fractional shortening and reduces aging-induced isovolumic relaxation time (IVRT) and cardiac hypertrophy in aged mice.

Reference

1. Kuppuswamy, D., Chinnakkannu, P., Reese, C., *et al.* The caveolin-1 scaffolding domain peptide reverses aging-associated deleterious changes in multiple organs. *J. Pharmacol. Exp. Ther.* **378**(1), 1-9 (2021).

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, 12/21/2022

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