

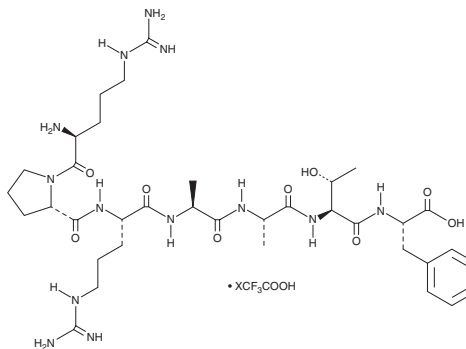
PRODUCT INFORMATION



Akt Substrate (trifluoroacetate salt)

Item No. 36759

Formal Name: L-arginyl-L-prolyl-L-arginyl-L-alanyl-L-alanyl-L-threonyl-L-phenylalanine, trifluoroacetate salt
Synonym: RPRAATF
MF: C₃₆H₅₉N₁₃O₉ • XCF₃COOH
FW: 817.9
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

Akt substrate (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Akt substrate (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Akt substrate (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

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 Akt substrate (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Akt substrate (trifluoroacetate salt) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Akt substrate is a peptide substrate for Akt, a serine/threonine kinase involved in various cellular processes such as proliferation, viability, angiogenesis, and glucose metabolism.¹ It has been used as a target substrate for the quantification of Akt activity.

Reference

1. Bozinovski, S., Cristiano, B.E., Marmy-Conus, N., *et al.* The synthetic peptide RPRAATF allows specific assay of Akt activity in cell lysates. *Anal. Biochem.* **305**(1), 32-39 (2002).

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