

PRODUCT INFORMATION



Biotin-SRC-1 (676-700) (trifluoroacetate salt)

Item No. 37511

Formal Name: N-[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]-L-cysteinyl-L-prolyl-L-seryl-L-seryl-L-histidyl-L-seryl-L-seryl-L-leucyl-L-threonyl-L- α -glutamyl-L-arginyl-L-histidyl-L-lysyl-L-isoleucyl-L-leucyl-L-histidyl-L-arginyl-L-leucyl-L-leucyl-L-glutamyl-L- α -glutamylglycyl-L-seryl-L-prolyl-L-serine, trifluoroacetate salt

Biotin—Cys—Pro—Ser—Ser—His—Ser—Ser—Leu—Thr—Glu—

Arg—His—Lys—Ile—Leu—His—Arg—Leu—Leu—Gln—

Glu—Gly—Ser—Pro—Ser—OH

• XCF₃COOH

Peptide Sequence: Biotin-CPSSHSLTERHKILHRLLEQEGSPS-OH
Synonym: Biotin-Steroid Receptor Coactivator 1 (676-700)

MF: C₁₂₈H₂₀₉N₄₁O₄₀S₂ • XCF₃COOH

FW: 3,026.4

Purity: ≥98%

Supplied as: A crystalline 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

Biotin-SRC-1 (676-700) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the biotin-SRC-1 (676-700) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Biotin-SRC-1 (676-700) is a biotinylated and truncated form of steroid receptor coactivator 1 (SRC-1). It has been used in affinity-based reporter assays, including time-resolved FRET (TR-FRET) and amplified luminescent proximity homogenous assay (ALPHASCREEN™).¹

Reference

1. Yang, F., Zhang, R., Ni, D., *et al.* Discovery of betulinaldehyde as a natural ROR γ t agonist. *Fitoterapia* 137, 104200 (2019).

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, 10/28/2024

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