

PRODUCT INFORMATION



BRC4wt (trifluoroacetate salt)

Item No. 38848

Formal Name: (S)-2-((5S,8S,11S,14S,17S,23S,26S)-14-((1H-imidazol-5-yl)methyl)-17-benzyl-11-((R)-1-hydroxyethyl)-5-(hydroxymethyl)-23,26-diisobutyl-8-methyl-4,7,10,13,16,19,22,25,28-nonaaxo-3,6,9,12,15,18,21,24,27-nonaazanacosanamido)-6-amino-N-((5S,8S,11S,14S,17S,20S)-1,24-diamino-14-(4-aminobutyl)-11-((S)-sec-butyl)-5-carbamoyl-17-isopropyl-8-methyl-7,10,13,16,19-pentaoxo-6,9,12,15,18-pentaazatetracosan-20-yl)hexanamide, trifluoroacetate salt

Ac—Leu—Leu—Gly—Phe—His—Thr—Ala—Ser—Gly—Lys—
Lys—Val—Lys—Ile—Ala—Lys—NH₂
• XCF₃COOH

Peptide Sequence: Ac-LLGFHTASGKKVKIAK-NH₂

MF: C₈₁H₁₃₉N₂₃O₁₉ • XCF₃COOH

FW: 1,739.1

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

BRC4wt (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the BRC4wt (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

BRC4wt is an acetylated peptide derived from the BRC4 repeat in human BRCA2 (1521-1536) and an inhibitor of the protein-protein interaction between BRCA2 and RAD51.¹ When coupled to the cationic cell-penetrating peptide (Arg)₉ (Item No. 36714), BRC4wt shortens DNA replication tract lengths and reduces homologous repair frequency of DNA damage induced by the DNA topoisomerase I inhibitor camptothecin (Item No. 11694) *in vitro*, as well as potentiates cell death induced by the poly(ADP-ribose) polymerase (PARP) inhibitor olaparib (Item No. 10621) in HeLa human cervical cancer and U2OS human osteosarcoma cells but not in non-cancerous hTERT RPE-1, MRC-5, or MCF-10A cells.

Reference

1. Trenner, A., Godau, J., and Sartori, A.A. A short BRCA2-derived cell-penetrating peptide targets RAD51 function and confers hypersensitivity toward PARP inhibition. *Mol. Cancer Ther.* **17(7)**, 1392-1404 (2018).

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, 07/28/2023

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