

PRODUCT INFORMATION

PD-1/PD-L1 Inhibitor 3 (trifluoroacetate salt)

Item No. 36719

Formal Name: cyclic (1→14)-thioether N-(2-mercaptoacetyl)-L-phenylalanyl-N-methyl-L-alanyl-L-asparaginyll-L-prolyl-L-histidyl-L-leucyl-N-methylglycyl-L-tryptophyl-L-seryl-L-tryptophyl-N-methyl-L-norleucyl-N-methyl-L-norleucyl-L-arginyl-L-cysteinyl-glycinamide

Synonym: Programmed Cell Death 1/Programmed Cell Death-Ligand 1 Inhibitor 3

MF: C₈₉H₁₂₆N₂₄O₁₈S • XCF₃COOH

FW: 1,852.2

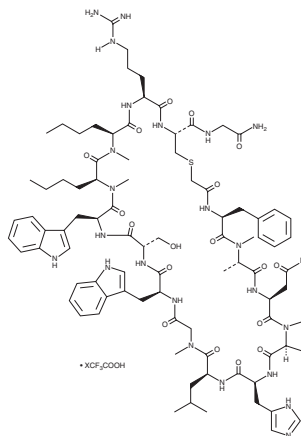
Purity: ≥95%

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

PD-1/PD-L1 inhibitor 3 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PD-1/PD-L1 inhibitor 3 (trifluoroacetate salt) in 30% acetonitrile/water. We do not recommend storing the aqueous solution for more than one day.

Description

PD-1/PD-L1 inhibitor 3 is a macrocyclic peptide inhibitor of the protein-protein interaction between programmed cell death protein 1 (PD-1) and PD-1 ligand (PD-L1).¹ It inhibits the PD-1/PD-L1 interaction in a time-resolved FRET (TR-FRET) assay (IC₅₀ = 9 nM).

Reference

1. Miller, M.M., Mapelli, C., Allen, P., *et al.* Macrocyclic inhibitors of the PD-1/PD-L1 and CD80(B7-1)/PD-L1 protein/protein interactions. *Bristol-Myers Squibb, Company US20140294898A1* (2017).

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