

PRODUCT INFORMATION

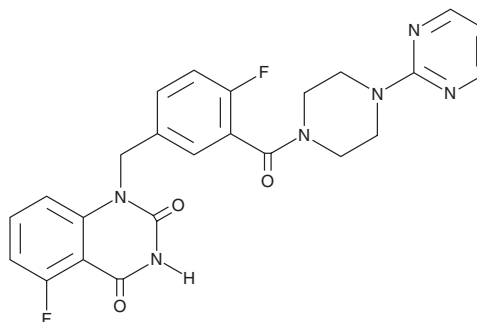


Senaparib

Item No. 37722

CAS Registry No.: 1401682-78-7
Formal Name: 5-fluoro-1-[[4-fluoro-3-[[4-(2-pyrimidinyl)-1-piperazinyl]carbonyl]phenyl]methyl]-2,4(1H,3H)-quinazolinedione

Synonym: IMP-4297
MF: C₂₄H₂₀F₂N₆O₃
FW: 478.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 244 nm
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

Senaparib is supplied as a solid. A stock solution may be made by dissolving the senaparib in the solvent of choice, which should be purged with an inert gas. Senaparib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of senaparib in ethanol is approximately 0.1 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

Senaparib is an inhibitor of poly(ADP-ribose) polymerase 1 (PARP1) and PARP2.¹

Reference

1. Meng, X., Lin, X., Jiang, R., *et al.* Effect of food on the pharmacokinetics of senaparib (IMP4297) in healthy Chinese subjects. *Clin. Drug Investig.* **42(11)**, 1009-1016 (2022).

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, 06/12/2023

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