

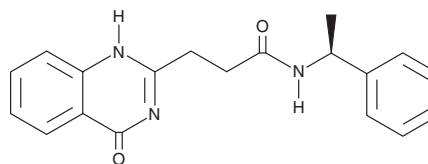
PRODUCT INFORMATION



ME0328

Item No. 18488

CAS Registry No.: 1445251-22-8
Formal Name: 3,4-dihydro-4-oxo-N-[(1S)-1-phenylethyl]-2-quinazolinepropanamide
MF: C₁₉H₁₉N₃O₂
FW: 321.4
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 263 nm
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

ME0328 is supplied as a crystalline solid. A stock solution may be made by dissolving the ME0328 in the solvent of choice, which should be purged with an inert gas. ME0328 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ME0328 in these solvents is approximately 0.2, 17, and 12 mg/ml, respectively.

ME0328 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ME0328 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ME0328 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ME0328 is an inhibitor of PARP3 (IC₅₀ = 0.89 μM), which is a regulator of DNA repair and mitotic progression.¹ It displays selectivity for PARP3 over 12 other PARP family members. ME0328 is cell permeable and elicits PARP3-specific effects at submicromolar concentrations.¹

Reference

1. Lindgren, A.E., Karlberg, T., Thorsell, A.G., *et al.* PARP inhibitor with selectivity toward ADP-ribosyltransferase ARTD3/PARP3. *ACS Chem. Biol.* **8**(8), 1698-1703 (2013).

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, 10/28/2022

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