

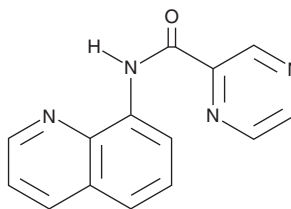
PRODUCT INFORMATION



QN523

Item No. 37123

CAS Registry No.: 878581-60-3
Formal Name: N-8-quinolinyl-2-pyrazinecarboxamide
MF: $C_{14}H_{10}N_4O$
FW: 250.3
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 240 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

QN523 is supplied as a crystalline solid. A stock solution may be made by dissolving the QN523 in the solvent of choice, which should be purged with an inert gas. QN523 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of QN523 in these solvents is approximately 2 and 5 mg/ml, respectively.

Description

QN523 is an anticancer agent.¹ It is cytotoxic to MiaPaCa-2, PANC-1, and BxPC-3 pancreatic cancer cells (IC_{50} s = 0.11, 0.5, and 3.3 μM , respectively). QN523 (0.5 μM) induces cell cycle arrest at the S phase in MiaPaCa-2 cells. It increases levels of the endoplasmic reticulum (ER) stress markers glucose-regulated protein 78 kDa (GRP78), DNA damage inducible transcript 3 (DDIT3), activating transcription factor 3 (ATF3), and growth differentiation factor 15 (GDF15), as well as the autophagy markers WD repeat domain phosphoinositide-interacting protein 1 (WIPI1), GABA_A receptor-associated protein-like 1 (GABARAPL1), and LC3-II in MiaPaCa-2 cells in a concentration-dependent manner. QN523 reduces tumor volume in a MiaPaCa-2 mouse xenograft model.

Reference

1. Kuang, Y., Ye, N., Kyani, A., *et al.* Induction of genes implicated in stress response and autophagy by a novel quinolin-8-yl-nicotinamide QN523 in pancreatic cancer. *J. Med. Chem.* **65**(8), 6133-6156 (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, 11/13/2025

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