

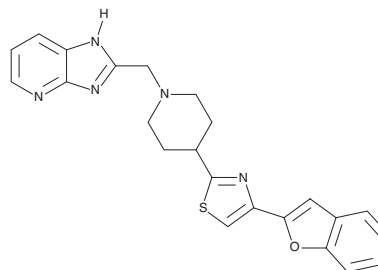
PRODUCT INFORMATION



BAY-179

Item No. 36657

CAS Registry No.: 2764880-87-5
Formal Name: 2-[[4-[4-(2-benzofuranyl)-2-thiazolyl]-1-piperidinyl]methyl]-3H-imidazo[4,5-b]pyridine
MF: C₂₃H₂₁N₅OS
FW: 415.5
Purity: ≥98%
UV/Vis.: λ_{max}: 286, 310 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

BAY-179 is supplied as a solid. A stock solution may be made by dissolving the BAY-179 in the solvent of choice, which should be purged with an inert gas. BAY-179 is slightly soluble in DMSO and dimethyl formamide.

Description

BAY-179 is an inhibitor of mitochondrial complex I, also known as NADH dehydrogenase (IC₅₀s = 79, 38, 27, and 47 nM for the human, mouse, rat, and dog enzymes, respectively).¹ See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Mowat, J., Ehrmann, A.H.M., Christian, S., *et al.* Identification of the highly active, species cross-reactive complex I inhibitor BAY-179. *ACS Med. Chem. Lett.* **13**(3), 348-357 (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, 05/08/2024

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