

PRODUCT INFORMATION

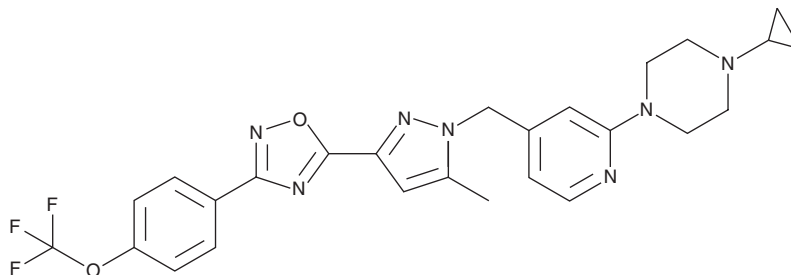


BAY 87-2243

Item No. 28626

CAS Registry No.: 1227158-85-1
Formal Name: 1-cyclopropyl-4-[4-[[5-methyl-3-[3-[4-(trifluoromethoxy)phenyl]-1,2,4-oxadiazol-5-yl]-1H-pyrazol-1-yl]methyl]-2-pyridinyl]-piperazine

MF: C₂₆H₂₆F₃N₇O₂
FW: 525.5
Purity: ≥98%
UV/Vis.: λ_{max}: 250 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

BAY 87-2243 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY 87-2243 in the solvent of choice, which should be purged with an inert gas. BAY 87-2243 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BAY 87-2243 in ethanol and DMSO is approximately 5 mg/ml and approximately 10 mg/ml in DMF.

BAY 87-2243 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAY 87-2243 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BAY 87-2243 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BAY 87-2243 is an inhibitor of mitochondrial complex I, also known as NADH:ubiquinone oxidoreductase, and an inducer of ferroptosis.^{1,2} It decreases ATP-dependent luciferase reporter activity and ATP levels in H1299luc cells, effects that can be reversed by addition of the mitochondrial complex II substrate succinate or expression of the *S. cerevisiae* complex I ortholog Ndi1 NADH dehydrogenase.¹ BAY 87-2243 induces production of reactive oxygen species (ROS), increases α-tocopherol-sensitive lipid peroxidation, and decreases glutathione (GSH) levels in SK-MEL-28 and G361 melanoma cells, effects that can be partially reversed by the ferroptosis inhibitor ferrostatin-1 (Item No. 17729) and overexpression of GPX4 or potentiated by GPX4 knockdown.² *In vivo*, BAY 87-2243 (0.5-4 mg/kg) reduces tumor weight, hypoxia-inducible factor-1α (HIF-1α) levels, and HIF-1α target gene expression in an H460 mouse xenograft model.¹ It also reduces tumor growth in the MEXF 276 and MEXF 1732 melanoma patient-derived xenograft (PDX) mouse models.³

References

1. Ellinghaus, P., Heisler, I., Unterschemmann, K., *et al.* BAY 87-2243, a highly potent and selective inhibitor of hypoxia-induced gene activation has antitumor activities by inhibition of mitochondrial complex I. *Cancer Med.* **2**(5), 611-624 (2013).
2. Basit, F., van Oppen, L.M., Schöckel, L., *et al.* Mitochondrial complex I inhibition triggers a mitophagy-dependent ROS increase leading to necroptosis and ferroptosis in melanoma cells. *Cell Death Dis.* **8**(3), e2716 (2017).
3. Schöckel, L., Glasauer, A., Basit, F., *et al.* Targeting mitochondrial complex I using BAY 87-2243 reduces melanoma tumor growth. *Cancer Metab.* **3**:11 (2015).

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, 10/27/2022

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