

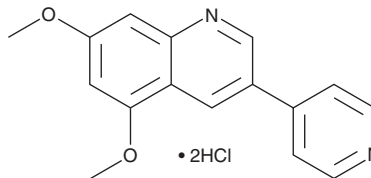
PRODUCT INFORMATION



DMPQ (hydrochloride)

Item No. 17551

CAS Registry No.: 1123491-15-5
Formal Name: 5,7-dimethoxy-3-(4-pyridinyl)-quinoline, dihydrochloride
MF: C₁₆H₁₄N₂O₂ • 2HCl
FW: 339.2
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 284, 374 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

DMPQ (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of DMPQ (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of DMPQ (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

DMPQ is a potent inhibitor of human platelet-derived growth factor receptor β (PDGFRβ, IC₅₀ = 80 nM).¹ It displays greater than 100-fold selectivity for PDGFRβ over other tyrosine and serine/threonine kinases.¹ DMPQ effectively blocks the PDGF-mediated increase in the protein, Survival Motor Neuron, in fibroblasts from patients with spinal muscular atrophy.²

References

1. Dolle, R.E., Dunn, J.A., Bobko, M., *et al.* 5,7-Dimethoxy-3-(4-pyridinyl)quinoline is a potent and selective inhibitor of human vascular β-type platelet-derived growth factor receptor tyrosine kinase. *J. Med. Chem.* **37**, 2627-2629 (1994).
2. Makhortova, N.R., Hayhurst, M., Cerqueira, A., *et al.* A screen for regulators of survival of motor neuron protein levels. *Nat. Chem. Biol.* **7**(8), 544-552 (2011).

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, 11/01/2022

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