

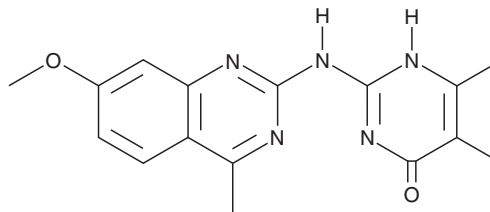
PRODUCT INFORMATION



Madrasin

Item No. 20451

CAS Registry No.: 374913-63-0
Formal Name: 2-[(7-methoxy-4-methyl-2-quinazolinyl) amino]-5,6-dimethyl-4(3H)-pyrimidinone
MF: C₁₆H₁₇N₅O₂
FW: 311.3
Purity: ≥95%
UV/Vis.: λ_{max}: 217, 253, 306, 330 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

Madrasin is supplied as a crystalline solid. A stock solution may be made by dissolving the madrasin in the solvent of choice. Madrasin is soluble in the organic solvent dimethyl formamide (DMF), which should be purged with an inert gas, at a concentration of approximately 1 mg/ml.

Madrasin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, madrasin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Madrasin has a solubility of approximately 0.025 mg/ml in a 1:40 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Madrasin is a cell-permeant inhibitor of pre-mRNA splicing.¹ It interferes with the early stages of spliceosome assembly and stalls spliceosome assembly at the A complex.¹ Madrasin induces cell cycle arrest, promotes a specific reorganization of subnuclear protein localization, and modulates splicing of multiple pre-RNA species in both HeLa and HEK293 cells when used at lower concentrations.¹

Reference

1. Pawellek, A., McElroy, S., Samatov, T., *et al.* Identification of small molecule inhibitors of pre-mRNA splicing. *J. Biol. Chem.* **289**(50), 34683-34698 (2014).

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, 11/14/2022

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