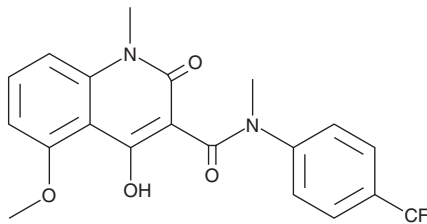


# PRODUCT INFORMATION



## Tasquinimod Item No. 17692

**CAS Registry No.:** 254964-60-8  
**Formal Name:** 1,2-dihydro-4-hydroxy-5-methoxy-N,1-dimethyl-2-oxo-N-[4-(trifluoromethyl)phenyl]-3-quinolinecarboxamide  
**Synonym:** ABR-215050  
**MF:** C<sub>20</sub>H<sub>17</sub>F<sub>3</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 406.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 232, 303 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

Tasquinimod is supplied as a crystalline solid. A stock solution may be made by dissolving the tasquinimod in the solvent of choice. Tasquinimod is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of tasquinimod in these solvents is approximately 2, 10, and 20 mg/ml, respectively.

Tasquinimod is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Tasquinimod is an orally-active quinoline-3-carboxamide derived from roquinimex, an immunomodulatory quinolone with applications in some cancers and autoimmune diseases.<sup>1</sup> Tasquinimod inhibits tumor angiogenesis and supplements radiation or chemotherapy in animal models of prostate cancer.<sup>1,2</sup> While its precise mechanism of action remains unclear, tasquinimod has been reported to alter signaling through S100A9, thrombospondin-1, HIF-1α, androgen receptor, VEGF, and HDAC3/4.<sup>1,2</sup>

### References

1. Isaacs, J.T. The long and winding road for the development of tasquinimod as an oral second-generation quinoline-3-carboxamide antiangiogenic drug for the treatment of prostate cancer. *Expert Opin. Investig. Drugs* **19**(10), 1235-1243 (2010).
2. Mehta, A.R. and Armstrong, A.J. Tasquinimod in the treatment of castrate-resistant prostate cancer - current status and future prospects. *Ther. Adv. Urol.* **8**(1), 9-18 (2016).

#### 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, 12/02/2022

#### 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](http://WWW.CAYMANCHEM.COM)