

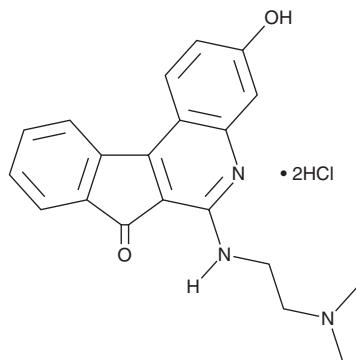
# PRODUCT INFORMATION



## TAS 103 (hydrochloride)

Item No. 29869

**CAS Registry No.:** 174634-09-4  
**Formal Name:** 6-[[2-(dimethylamino)ethyl]amino]-3-hydroxy-7H-indeno[2,1-c]quinolin-7-one, dihydrochloride  
**MF:**  $C_{20}H_{19}N_3O_2 \cdot 2HCl$   
**FW:** 406.3  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 229, 244, 262, 292 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

TAS 103 (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of TAS 103 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TAS 103 (hydrochloride) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

TAS 103 is an inhibitor of DNA topoisomerase I and II ( $IC_{50}s = 2$  and  $6.5 \mu M$ , respectively).<sup>1</sup> It stabilizes topoisomerase I-DNA and topoisomerase II-DNA cleavable complexes in KB cells when used at a concentration of  $3 \mu M$ . It is cytotoxic to P388 and KB cells ( $IC_{50}s = 1.1$  and  $9.6 nM$ , respectively). TAS 103 (15, 26, and  $45 mg/kg$  per day) reduces pulmonary metastasis in murine B16-BL6 melanoma and UV-2237M fibrosarcoma models of lung metastasis. It also reduces tumor growth in a variety of human cancer mouse xenograft models.

### Reference

1. Utsugi, T., Aoyagi, K., Asao, T., *et al.* Antitumor activity of a novel quinoline derivative, TAS-103, with inhibitory effects on topoisomerases I and II. *Jpn. J. Cancer Res.* **88**(10), 992-1002 (1997).

#### 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, 12/12/2022

#### 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)