

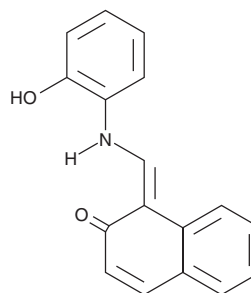
PRODUCT INFORMATION



HAMNO

Item No. 27631

CAS Registry No.: 138736-73-9
Formal Name: 1-[[[(2-hydroxyphenyl)amino]methylene]-2(1H)-naphthalenone
MF: C₁₇H₁₃NO₂
FW: 263.3
Purity: ≥95%
UV/Vis.: λ_{max}: 211, 235, 263, 321, 448, 472 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

HAMNO is supplied as a crystalline solid. A stock solution may be made by dissolving the HAMNO in the solvent of choice, which should be purged with an inert gas. HAMNO is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of HAMNO in these solvents is approximately 30 mg/ml.

HAMNO is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HAMNO should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. HAMNO has a solubility of approximately 0.25 mg/ml in a 1:30 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HAMNO is a small molecule inhibitor of replication protein A (RPA).¹ It selectively binds to and inhibits DNA binding domain F (DBD-F) on RPA70 when used at concentrations up to 100 μM. HAMNO inhibits etoposide-induced autophosphorylation of ATR kinase and subsequent phosphorylation of RPA32 in UMSCC38 cells when used at concentrations of 10 and 20 μM. HAMNO (1 mg/kg), in combination with etoposide (Item No. 12092), reduces tumor growth in UMSCC11B and UMSCC38 mouse xenograft models.

Reference

1. Glanzer, J.G., Liu, S., Wang, L., *et al.* RPA inhibition increases replication stress and suppresses tumor growth. *Cancer Res.* **74(18)**, 5165-5172 (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, 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