

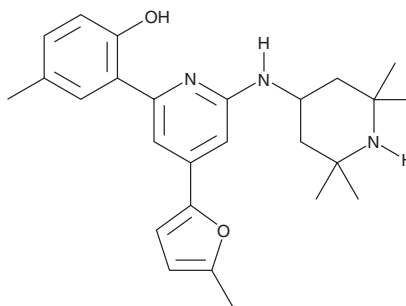
PRODUCT INFORMATION



BRITE338733

Item No. 37302

CAS Registry No.: 503105-88-2
Formal Name: 2-[4-(5-ethyl-2-furanyl)-6-[(2,2,6,6-tetramethyl-4-piperidiny)amino]-2-pyridinyl]-4-methyl-phenol
MF: C₂₇H₃₅N₃O₂
FW: 433.6
Purity: ≥95%
UV/Vis.: λ_{max}: 211, 290 nm
Supplied as: A 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

BRITE338733 is supplied as a solid. A stock solution may be made by dissolving the BRITE338733 in the solvent of choice, which should be purged with an inert gas. BRITE338733 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRITE338733 in ethanol and DMF is approximately 1 mg/ml and approximately 3 mg/ml in DMSO.

Description

BRITE338733 is an inhibitor of *E. coli* recombinase A (RecA; IC₅₀ = 4.7 μM for ATP hydrolysis in a cell-free assay), an enzyme involved in DNA repair.¹

Reference

1. Sexton, J.Z., Wigle, T.J., He, Q., *et al.* Novel inhibitors of *E. coli* RecA ATPase activity. *Curr. Chem. Genomics* 4, 34-42 (2010).

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, 04/10/2026

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