

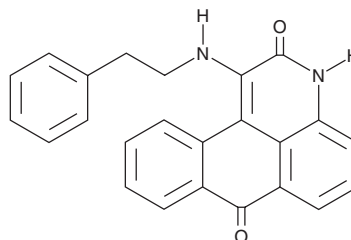
PRODUCT INFORMATION



BRD7389

Item No. 20214

CAS Registry No.: 376382-11-5
Formal Name: 1-[(2-phenylethyl)amino]-3H-naphtho[1,2,3-de]quinoline-2,7-dione
MF: C₂₄H₁₈N₂O₂
FW: 366.4
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 239, 343, 444 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

BRD7389 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD7389 in the solvent of choice, which should be purged with an inert gas. BRD7389 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BRD7389 in these solvents is approximately 12.5 and 20 mg/ml, respectively. BRD7389 is slightly soluble in ethanol.

BRD7389 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD7389 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BRD7389 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD7389 is an inhibitor of ribosomal S6 kinases (RSKs; IC₅₀s = 1.5, 2.4, and 1.2 μM for RSK1, RSK2, and RSK3, respectively).¹ Through this action, BRD7389 causes mouse αTC1 pancreatic α cells to adopt several morphological and gene expression features of β cells, including increased insulin expression.¹ BRD7389 also enhances glucose-stimulated insulin and glucagon secretion by pancreatic islet cells isolated from human donors with a low body mass index.¹

Reference

1. Fomina-Yadlin, D., Kubicek, S., Walpita, D., *et al.* Small-molecule inducers of insulin expression in pancreatic α-cells. *Proc. Natl. Acad. Sci. USA* **107**(34), 15099-15104 (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, 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