

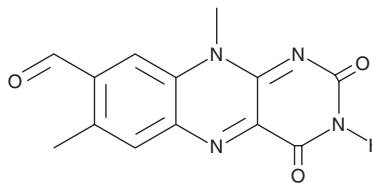
PRODUCT INFORMATION



Ro 08-2750

Item No. 28772

CAS Registry No.: 37854-59-4
Formal Name: 2,3,4,10-tetrahydro-7,10-dimethyl-2,4-dioxo-benzo[g]pteridine-8-carboxaldehyde
MF: C₁₃H₁₀N₄O₃
FW: 270.2
Purity: ≥95%
UV/Vis.: λ_{max}: 248, 288, 335, 460 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

Ro 08-2750 is supplied as a solid. A stock solution may be made by dissolving the Ro 08-2750 in the solvent of choice, which should be purged with an inert gas. Ro 08-2750 is soluble in the organic solvent DMSO.

Description

Ro 08-2750 is a p75 neurotrophin receptor (p75^{NTR}) ligand that completely inhibits nerve growth factor (NGF) binding to p75^{NTR} in PC12 cells when used at a concentration of 100 μM.¹ It also binds to NGF dimers (K_d = ~1 μM) to induce a conformational change that prevents NGF binding to p75^{NTR} or tropomyosin-related kinase A (TrkA). Ro 08-2750 inhibits NGF-induced death of SK-N-MC cells and inhibits autophosphorylation of TrkA in PC12 cells. It reverses NGF-induced decreases in β-catenin mRNA expression in, and migration of, SKOV3 and OVCAR-3 ovarian cancer cells.²

References

1. Niederhauser, O., Mangold, M., Schubeneil, R., *et al.* NGF ligand alters NGF signaling via p75^{NTR} and trkA. *J. Neurosci. Res.* **61(3)**, 263-272 (2000).
2. Li, B., Cai, S., Zhao, Y., *et al.* Nerve growth factor modulates the tumor cells migration in ovarian cancer through the WNT/β-catenin pathway. *Oncotarget* **7(49)**, 81026-81048 (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, 12/02/2019

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