

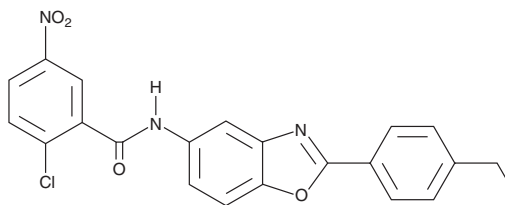
PRODUCT INFORMATION



BAY-4931

Item No. 40668

CAS Registry No.: 423150-91-8
Formal Name: 2-chloro-N-[2-(4-ethylphenyl)-5-benzoxazolyl]-5-nitro-benzamide
MF: $C_{22}H_{16}ClN_3O_4$
FW: 421.8
Purity: $\geq 98\%$
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

BAY-4931 is supplied as a solid. A stock solution may be made by dissolving the BAY-4931 in the solvent of choice, which should be purged with an inert gas. BAY-4931 is soluble (≥ 10 mg/ml) in DMSO.

Description

BAY-4931 is a PPAR γ inverse agonist.¹ It acts as an inverse agonist in a co-repressor reporter assay with a peptide from nuclear receptor co-repressor 2 (NCOR2; $EC_{50} = 5.8$ nM). BAY-4931 selectively induces PPAR γ -mediated gene transcription ($IC_{50} = 0.4$ nM for the human receptor) over PPAR α - or PPAR δ -mediated gene transcription in reporter assays using CHO cells (IC_{50} s = >50 μM for both for the human receptors). BAY-4931 inhibits the proliferation of UM-UC-9 human bladder cancer cells that endogenously contain a focal amplification of *PPARG*, the gene encoding PPAR γ ($IC_{50} = 3.4$ nM).

Reference

- Orsi, D.L., Pook, E., Bräuer, N., *et al.* Discovery and structure-based design of potent covalent PPAR γ inverse-agonists BAY-4931 and BAY-0069. *J. Med. Chem.* **65**(21), 14843-14863 (2022).

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, 09/05/2024

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