PNU-74654

Catenin blocker

Anti-tumor agent that disrupts the canonical Wnt/ β -catenin pathway *via* blockage of the protein-protein interaction between β -catenin and TCF.

Citations: 3

View Online »

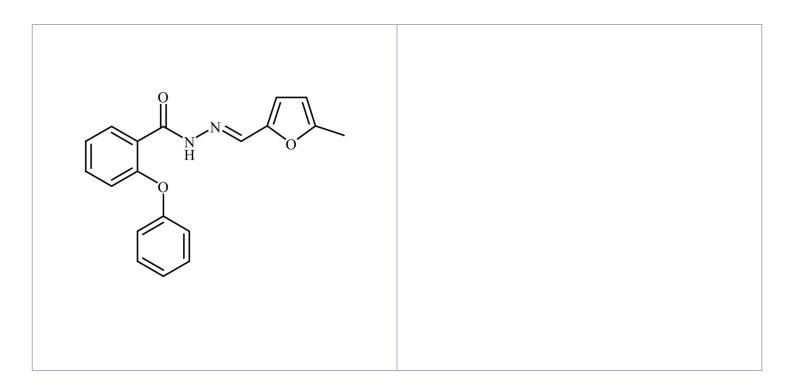
Ordering Information

Order Online »

BML-WN101-0005	5mg
BML-WN101-0025	25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

6 months after receipt when stored at -20°C.

Handling Solutions can be stored at -20°C for up to 3 months.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-Phenoxybenzoic acid [(5-methyl-2-

furanyl)methylene]hydrazide

Appearance Off-white solid.

CAS 113906-27-7

Couple Target Catenin

Couple Type Blocker

Formula $C_{19}H_{16}N_2O_3$

MW 320.3

Purity ≥98% (HPLC)

Soluble in DMSO (25 mg/ml) and in 100% ethanol (8

mg/ml).

Last modified: May 29, 2024

