

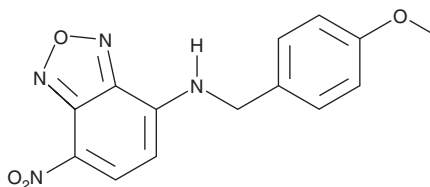
PRODUCT INFORMATION



7-(*p*-Methoxybenzylamino)-4-nitrobenz-2,1,3-oxadiazole

Item No. 39668

CAS Registry No.: 33984-50-8
Formal Name: N-[(4-methoxyphenyl)methyl]-7-nitro-2,1,3-benzoxadiazol-4-amine
Synonyms: NSC 240872, MBD
MF: C₁₄H₁₂N₄O₄
FW: 300.3
Purity: ≥98%
Ex./Em. Max: 430/530 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

7-(*p*-Methoxybenzylamino)-4-nitrobenz-2,1,3-oxadiazole (MBD) is supplied as a solid. A stock solution may be made by dissolving the MBD in the solvent of choice, which should be purged with an inert gas. MBD is soluble in methanol and DMSO.

Description

MBD is a fluorescent probe for hydrophobic regions of proteins.^{1,2} Upon binding to a hydrophobic region of a protein, MBD displays an emission of 530 nm upon excitation at 430 nm.¹

References

1. Kenner, R.A. and Aboderin, A.A. A new fluorescent probe for protein and nucleoprotein conformation. Binding of 7-(*p*-methoxybenzylamino)-4-nitrobenzoxadiazole to bovine trypsinogen and bacterial ribosomes. *Biochemistry* **10**(24), 4433-4440 (1971).
2. Kasahara, Y., Takahashi, K., Nagasawa, S., *et al.* Formation of a conformationally changed C1r, a subcomponent of the first component of human complement, as an intermediate of its autoactivation reaction. *FEBS Lett.* **141**(1), 128-131 (1982).

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, 02/19/2024

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