

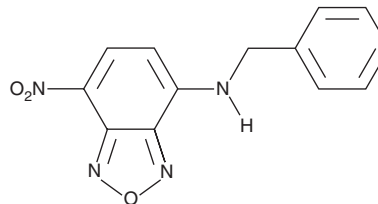
PRODUCT INFORMATION



BBD

Item No. 42897

CAS Registry No.: 18378-20-6
Formal Name: 7-nitro-N-(phenylmethyl)-2,1,3-benzoxadiazol-4-amine
Synonyms: Benzylamino-NBD, 4-Benzylamino-7-nitrobenzofurazan, NSC 240867
MF: $C_{13}H_{10}N_4O_3$
FW: 270.2
Purity: $\geq 95\%$
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

BBD is supplied as a solid. A stock solution may be made by dissolving the BBD in the solvent of choice, which should be purged with an inert gas. BBD is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

BBD is a fluorescent probe for the detection of hydrophobicity.¹⁻³ It exhibits absorption and emission maxima that shift to shorter wavelengths in response to increasing concentrations of ethanol in water.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. Stevenson, S.A. and Blanchard, G.J. Dynamics of 4-benzylamino-7-nitrobenzofurazan in the 1-propanol/water binary solvent system. Evidence for composition-dependent solvent organization. *J. Phys. Chem. A* **110**(10), 3426-3431 (2006).
3. Matsushita, Y., Takahashi, M., and Moriguchi, I. Binding of fluorescent 7-amino-4-nitrobenzoxadiazole derivatives to bovine serum albumin. *Chem. Pharm. Bull. (Tokyo)* **34**(1), 333-339 (1986).

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, 02/20/2025

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