

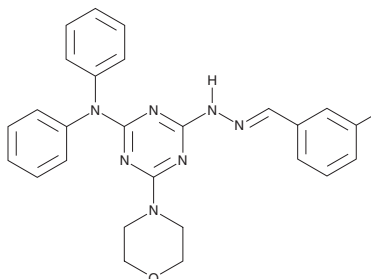
PRODUCT INFORMATION



Vacuolin-1

Item No. 20425

CAS Registry No.: 351986-85-1
Formal Name: 3-iodo-benzaldehyde,2-[4-(diphenylamino)-6-(4-morpholinyl)-1,3,5-triazin-2-yl]hydrazone
MF: C₂₆H₂₄IN₇O
FW: 577.4
Purity: ≥95%
UV/Vis.: λ_{max}: 225, 242, 314 nm
Supplied as: A crystalline 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

Vacuolin-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the vacuolin-1 in the solvent of choice, which should be purged with an inert gas. Vacuolin-1 is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

Vacuolin-1 is a cell-permeable inhibitor of lysosomal exocytosis that prevents Ca²⁺-dependent fusion of lysosomes with the plasma membrane and the release of lysosomal content.¹ It does not affect other membrane-bound organelles (including enlargeosomal exocytosis), disturb the organization of the actin or tubulin cytoskeletal network, nor interfere with trafficking along the endocytic or secretory pathways.¹

Reference

1. Cerny, J., Feng, Y., Yu, A., *et al.* The small chemical vacuolin-1 inhibits Ca²⁺-dependent lysosomal exocytosis but not cell resealing. *EMBO Rep.* 5(9), 883-888 (2004).

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, 07/06/2023

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