

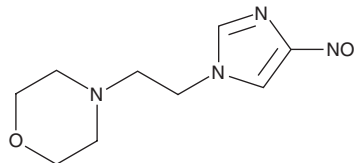
PRODUCT INFORMATION



RGW 611

Item No. 30067

CAS Registry No.: 6497-78-5
Formal Name: 4-[2-(4-nitro-1H-imidazol-1-yl)ethyl]-morpholine
MF: C₉H₁₄N₄O₃
FW: 226.2
Purity: ≥98%
UV/Vis.: λ_{max}: 298 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

RGW 611 is supplied as a crystalline solid. A stock solution may be made by dissolving the RGW 611 in the solvent of choice, which should be purged with an inert gas. RGW 611 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of RGW 611 in these solvents is approximately 1 and 5 mg/ml, respectively.

Description

RGW 611 is a nitroimidazole.¹ It enhances radiation-induced cell death of hypoxic V79-379A cells.

Reference

1. Adams, G.E., Clarke, E.D., Flockhart, I.R., *et al.* Structure-activity relationships in the development of hypoxic cell radiosensitizers. I. Sensitization efficiency. *Int. J. Radiat. Biol. Relat. Stud. Phys. Chem. Med.* **35**(2), 133-150 (1979).

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, 05/12/2020

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