

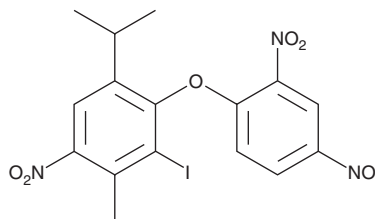
PRODUCT INFORMATION



DNP-INT

Item No. 9003061

CAS Registry No.: 69311-70-2
Formal Name: 2-(2,4-dinitrophenoxy)-3-iodo-4-methyl-1-(1-methylethyl)-5-nitro-benzene
MF: C₁₆H₁₄IN₃O₇
FW: 487.2
Purity: ≥98%
UV/Vis.: λ_{max}: 280 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

DNP-INT is supplied as a crystalline solid. A stock solution may be made by dissolving the DNP-INT in the solvent of choice. DNP-INT is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of DNP-INT in these solvents is approximately 5 and 10 mg/ml, respectively.

DNP-INT is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DNP-INT should first be dissolved in DMF and then diluted with the aqueous buffer of choice. DNP-INT has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DNP-INT is a quinone analog that inhibits electron transport in plants by competitively inhibiting plastoquinol oxidation by binding at the Q_o site of cytochrome *b₆f* (K_d = 1.4 nM).¹ It inhibits electron flow from water to NADP or methylviologen by 50 and 100% when used at concentrations of 0.5 or 5 μM, respectively.^{2,3}

References

1. Barbagallo, R.P., Finazzi, G., and Forti, G. Effects of inhibitors on the activity of the cytochrome *b₆f* complex: Evidence for the existence of two binding pockets in the lumenal site. *Biochemistry* **38**(39), 12814-12821 (1999).
2. Trebst, A., Wietoska, H., Draber, W., *et al.* The inhibition of photosynthetic electron flow in chloroplasts by the dinitrophenylether of bromo- or iodo-nitrothymol. *Zeitschrift für Naturforschung C* **33**(11-12), 919-927 (1978).
3. Malkin, R. Interaction of stigmatellin and DNP-INT with the Rieske iron-sulfur center of the chloroplast cytochrome *b₆f* complex. *FEBS Lett.* **208**(2), 317-320 (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, 10/17/2017

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