

DNP maleimide

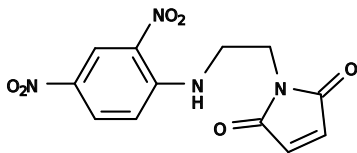
Catalog Number: 2026

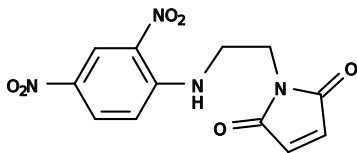
Unit Size: 25 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid yellow
Molecular Weight	306.23
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

A thiol-reactive building block for developing a probe that can be recognized by anti-DNP antibodies; Also an excellent thiol-reactive FRET quencher paired with Trp or Tyr.