

Product Information Sheet

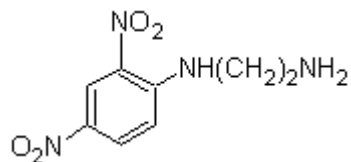
Ordering Information

Product Number: 2025
Product Name: DNP amine
Unit Size: 25 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 226.19

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: 350 nm

Emission Wavelength: NA

Application Notes

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