

## DNP-X acid, SE [6-(2,4-Dinitrophenyl)amino]hexanoic acid, succinimidyl ester] \*CAS 82321-04-8\*

Catalog Number: 2021

Unit Size: 25 mg

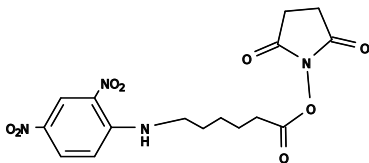
### Product Details

---

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

### Chemical Properties

---

Appearance	Solid yellow
Molecular Weight	394.34
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

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