

## DNP-X acid [6-(2,4-Dinitrophenyl)aminohexanoic acid] \*CAS 10466-72-5\*

Catalog Number: 2020

Unit Size: 100 mg

### Product Details

---

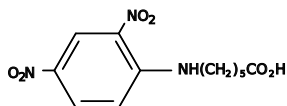
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	297.26
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.