

DNP-dPEG4-NHS Ester

A DNP moiety containing a PEG linker and an NHS ester terminal group which can be used for labeling biomolecules with a DNP hapten tag. The NHS ester will react with primary amines to form a stable amide bond while the hydrophilic PEG linker improves the solubility of the labeled bioconjugate.



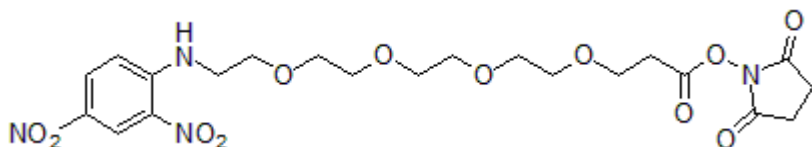
Product attributes

CAS number	858126-78-0
------------	-------------

Product Description

DNP is a proton ionophore that is often used in biochemistry as a hapten tag that can be detected using an anti-DNP antibody. DNP-dPEG4-NHS Ester is a DNP moiety containing a PEG linker and an NHS ester terminal group which can be used for labeling biomolecules. The NHS ester will react with primary amines to form a stable amide bond while the hydrophilic PEG linker improves the solubility of the labeled bioconjugate. DNP-dPEG4-NHS Ester may also be used as an amine-reactive FRET quencher when paired with tryptophan or tyramine.

- Yellow solid soluble in DMSO or DMF
- Supplied as 50 mg or 500 mg
- Store at -20°C and protect from light
- $C_{21}H_{28}N_4O_{12}$
- MW: 529
- [858126-78-0]



This datasheet was generated on June 7, 2026 at 04:37:17 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dnp-dpeg4-nhs-ester/>