## **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

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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<sub>21</sub>H<sub>28</sub>N<sub>4</sub>O<sub>12</sub>
- MW: 529
- [858126-78-0]

$$NO_2$$
 $NO_2$ 
 $NO_2$ 
 $NO_2$ 
 $NO_2$ 

This datasheet was generated on November 2, 2025 at 08:58:31 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/dnp-dpeg4-nhs-ester/">https://biotium.com/product/dnp-dpeg4-nhs-ester/</a>