

Product Information

DNP-dPEG4-NHS Ester

Catalog Number: 91062-50mg, 91062-500mg

Unit Size: 50 mg, 500 mg

Storage and Handling

Store at -20°C desiccated and protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

Stock solutions may be prepared in DMSO or DMF. Solutions can be aliquoted and stored with desiccant and protected from light at -20°C, for up to 12 months.

Molecular Information: $C_{21}H_{28}N_4O_{12}$

CAS number: 858126-78-0 Molecular Weight: 529 Color and Form: Yellow solid Solubility: DMSO or DMF

$$\begin{array}{c|c}
H \\
NO_2
\end{array}$$

Figure 1. DNP-dPEG4-NHS Ester

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.

Related Products

Cat. No.	Product
90082	DMSO, Anhydrous
92103 97500	CF® Dye SE/TFP Ester
92325- 92327	Mix-n-Stain™ DNP Antibody Labeling Kit
96019	DNP Tyramide
20865 20875	DNP Polyclonal Rabbit Antibody

Please visit our website at www.biotium.com for information on our life science research products, including rapid and efficient Mix-n-Stain™ Antibody Labeling Kits, master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.