

# Tide Quencher™ 7HWS Succinimidyl Ester [TQ7HWS Succinimidyl Ester]

Catalog Number: 2389

Unit Size: 1 mg

## Product Details

---

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

## Chemical Properties

---

Appearance	Solid
Molecular Weight	1391.84
Soluble In	DMSO

## Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

---

The non-fluorescent Tide Quencher™ 7HWS Succinimidyl Ester (TQ7HWS Succinimidyl Ester) is a next-generation, highly water-soluble quencher optimized for fluorescence-based applications, particularly in oligonucleotide and peptide labeling. As an advanced derivative of TQ7WS, it features significantly enhanced water solubility, enabling efficient labeling in aqueous environments with minimal organic solvent requirements. Like TQ7WS, TQ7HWS is designed to be a superior quencher for fluorophores such as Tide Fluor™ 7 (TF7), Alexa Fluor 750, and Cy7, offering stronger absorption properties, higher quenching efficiency, and excellent solubility. The increased hydrophilicity of TQ7HWS acid further improves its applicability in biological and biochemical assays requiring stringent aqueous conditions. With its exceptional quenching performance, complete absence of background fluorescence, and excellent solubility, TQ7HWS provides a reliable and efficient solution for demanding fluorescence assays.

Tide Quencher™ 7HWS Succinimidyl Ester (TQ7HWS Succinimidyl Ester) selectively and efficiently reacts with primary amines, including lysine side chains in peptides and aminosilane-coated surfaces, under mildly basic conditions (pH 7–9). This reaction forms stable covalent amide bonds, making TQ7HWS Succinimidyl Ester a valuable reagent for peptide labeling. Additionally, it is utilized for the post-synthetic labeling of amino-modified oligonucleotides.