

## Tide Quencher™ 5HWS Acid [TQ5HWS Acid]

Catalog Number: 2384

Unit Size: 5 mg

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	1078.17
Soluble In	DMSO

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

The non-fluorescent Tide Quencher™ 5HWS Acid (TQ5HWS acid) is a next-generation, highly water-soluble quencher optimized for fluorescence-based applications, particularly in oligonucleotide and peptide labeling. As an advanced derivative of TQ5WS, it features significantly enhanced water solubility, enabling efficient labeling in aqueous environments with minimal organic solvent requirements. Like TQ5WS, TQ5HWS is designed to be a superior quencher for fluorophores such as Tide Fluor™ 5 (TF5), Cy5, and Cy5.5, offering stronger absorption properties, higher quenching efficiency, and excellent solubility. The increased hydrophilicity of TQ5HWS 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, TQ5HWS provides a reliable and efficient solution for demanding fluorescence assays.

TQ5HWS acid is a non-reactive compound used as a reference standard or control in studies involving TQ5HWS-labeled conjugates. Additionally, it serves as a key precursor for the synthesis of activated esters and STP derivatives, facilitating further chemical modifications. In aqueous solutions, TQ5HWS acid can be conjugated to hydrazines, hydroxylamines, or amines through water-soluble carbodiimides, such as EDAC. This enables the attachment of the quencher to amino-containing biomolecules, including amine-modified oligonucleotides and peptides.