

Tide Quencher™ 2WS Azide [TQ2WS Azide]

Catalog Number: 2215

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	36 months upon receiving

Chemical Properties

Appearance	Solid dark red
Molecular Weight	1012.29
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Tide Quencher™ 2WS Azide is a highly water-soluble dark quencher engineered for efficient and specific conjugation to alkyne-modified biomolecules through azide-based click chemistry. It is ideal for labeling oligonucleotides, peptides, and other click-compatible molecules for use in molecular biology, FRET-based assays, and bioconjugation workflows.

TQ2WS Azide exhibits strong absorption and efficient quenching across a broad spectral range, effectively quenching dyes such as FAM, HEX, TET, JOE, TF2, and Rhodamine 6G. Its non-fluorescent profile minimizes background interference, supporting applications such as molecular probes and qPCR. Compared to standard TQ2, the enhanced water solubility of TQ2WS Azide is particularly beneficial for improving the solubility of hydrophobic targets during synthesis and labeling.