

# Tide Quencher™ 7.2 CPG [TQ7.2 CPG] \*1000 Å\*

Catalog Number: 2125

Unit Size: 100 mg

## Product Details

---

Storage Conditions                      Freeze (< -15 °C), Minimize light exposure

Expiration Date                          12 months upon receiving

## Chemical Properties

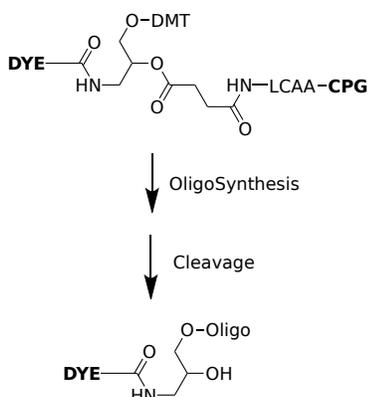
---

Appearance                                Solid green

Molecular Weight                        N/A

Soluble In                                  MeCN

Chemical Structure



## Spectral Properties

---

Excitation Wavelength                  N/A

Emission Wavelength                  N/A

## Applications

---

Tide Quencher™ 7.2 (TQ7.2) is a non-fluorescent molecule designed to efficiently quench the fluorescence of common NIR fluorophores such as Cy7, Alexa Fluor® 700, Alexa Fluor® 750, iFluor® 700, iFluor® 710, iFluor® 720 and iFluor® 750. It is an improved version of TQ7 and BHQ3. TQ7.2 is designed to be a superior quencher with (a). much stronger absorption, and (b). much higher quenching efficiency for NIR dyes. Tide Quencher™ 7.2 CPG is an excellent building block for preparing TQ7.2-labeled oligonucleotides. The oligo prepared from (TQ7.2 CPG) may be deprotected in 0.05M potassium carbonate in methanol for 4 hours at room temperature for 2 hours. Alternatively, the oligo may be deprotected in ammonium hydroxide at room temperature for 24-36 hours. It can be used in techniques such as polymerase chain reaction (PCR), real-time PCR, and DNA sequencing. In these applications, fluorescence signals are used to monitor the amplification or detection of specific DNA sequences. TQ7.2 quenches the fluorescent signal until a specific event (like DNA strand separation or primer extension) occurs, leading to an increase in fluorescence that can be detected and quantified.