

Texas Red™ Maleimide

Cat. No. SP-1505

Storage Store at -20 to -80 °C in the dark.

5' EndTag™ or 3' EndTag™ labeling of nucleic acids

Dissolve Texas Red Maleimide with 500 µl of anhydrous dimethyl sulfoxide (DMSO) and store at -20 °C to -80 °C in the dark.

Follow the labeling procedure included with the 5' EndTag or 3' EndTag systems.

Protein labeling procedure

1. Dissolve the protein to be labeled in 0.1 M phosphate buffer, 5 mM EDTA, pH 7, at a concentration of 5 mg/ml.
2. Dissolve a slight excess of the Texas Red Maleimide needed in dimethyl sulfoxide (DMSO) at 20 mg/ml.
3. Add 25 µl Texas Red Maleimide per ml of protein solution.
4. Incubate at room temperature for 3 hours with occasional stirring.
5. Separate the unreacted material from the protein by gel filtration or dialysis.

References

¹Deziel, M.R. et al. 1990. Biotin-conjugated reagents as site-specific probes of membrane protein structure: application to the study of the human erythrocyte hexose transporter. *Anal. Biochem.*

²Sambrook, J. et al. 1989. *Molecular Cloning: A Laboratory Manual*. Cold Spring Harbor Laboratory Press