

# Fluorescein Maleimide

**Cat. No.** SP-1502

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

## 3' EndTag™ labeling of nucleic acids

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

Follow the labeling procedure included with the 3' EndTag system.

## 5' EndTag™ labeling of nucleic acids

Dissolve Fluorescein Maleimide with 883 µl of anhydrous dimethyl sulfoxide (DMSO) and store at -20 to -80 °C in the dark.

Follow the labeling procedure included with the 5' EndTag system.

## Protein labeling procedure

1. Dissolve the protein to be labeled in 100 mM phosphate buffer, pH 7.0 at a concentration of 5 mg/ml.
2. Dissolve a slight excess of the amount needed of Fluorescein Maleimide in dimethyl Sulfoxide (DMSO) at a concentration of 20 mg/ml.
3. Add 25 µl Fluorescein 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

<sup>1</sup>Deziel, 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.*

## Selected reagents for the detection of the fluorescein label

Goat Anti-Fluorescein, Biotinylated	BA-0601	0.5 mg
-------------------------------------	---------	--------