

# Biotin (Long Arm) Maleimide

**Cat. No.** SP-1501

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

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

Dissolve Biotin (Long Arm) Maleimide in 312 µl of anhydrous dimethyl formamide (DMF) 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 100 mM phosphate buffer, pH 7.0 at a concentration of 5 mg/ml.
2. Dissolve a slight excess of the amount needed of Biotin (Long Arm) Maleimide in dimethyl formamide (DMF) at a concentration of 20 mg/ml.
3. Add 25 µl Biotin (Long Arm) 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. and Mau, M.M. 1990. Biotin-conjugated reagents as site-specific probes of membrane protein structure: application to the study of the human erythrocyte hexose transporter. *Anal. Biochem.* 190:297-303

## Selected reagents for the detection of the biotin label

Streptavidin, Alkaline Phosphatase	SA-5100	1 ml
Anti-Biotin Mouse Monoclonal	MB-9100	1 ml
Goat Anti-Biotin, Unconjugated	SP-3000	1 mg
Goat Anti-Biotin, Peroxidase	SP-3010	1 mg
Streptavidin, Peroxidase	SA-5004	1 mg
VECTASTAIN® ABC Kit (Standard)	PK-6100	1 kit