

---

## Product Information Sheet

### **Ordering Information**

Product Number: 11026  
Product Name: ReadiUse™ Preactivated HRP maleimide  
Unit Size: 300 ug  
Storage Conditions: Freeze (<-15 °C)  
Expiration Date: 12 months upon receiving

### **Chemical and Spectral Properties**

Appearance: Solid  
Molecular Weight: N/A  
Soluble In: N/A  
Excitation Wavelength: N/A  
Emission Wavelength: N/A

### **Application Notes**

ReadiUse™ Preactivated maleimide is a thiol-reactive HRP derivative. It can be used to readily label proteins (such as antibodies) or other biological molecules that have a thiol group. ReadiUse™ Preactivated HRP maleimide is robust and easy to use. It can be used to label antibodies with a simple mixing, i.e. the slightly acidic or neutral solution of an antibody (pH 6-7) can be directly mixed with the HRP maleimide, and the reaction mixture is shaken for 1-2 hours. In most cases, the resulted solution can be directly used for ELISA assays without further purification.