

ReadiUse™ Preactivated HRP NHS Ester (Cat# 11025, 300 ug)

Ordering Information:

Product Number: 11025 (300 ug)

Storage Conditions:

Keep at -20 °C
Avoid exposure to light

Introduction

ReadiUse™ Preactivated HRP NHS ester is the mono NHS ester of HRP. It can be used to readily label proteins (such as antibodies) or other biologicals molecules that have an amino group. ReadiUse™ Preactivated HRP NHS ester is robust and easy to use. It can be used to label antibodies (~ 100 ug) with a simple mixing, i.e. the basic solution of an antibody (pH 8.5-9.0) can be directly mixed with the HRP NHS ester (300 ug). In most cases, the resulted solution can be directly used for ELISA assays without further purification.

Assay Protocol:

Note 1: If the target protein solution contains other proteins (such as BSA), purify the protein solution or it will interfere with the labeling reaction.

Note 2: Small molecules such as sodium azide need to be removed.

Note 3: Tris buffer and other amine compounds will also interfere with the labeling reaction.

1. Antibody (IgG) ~100 µg in 20 µL~50µL of 25mM Sodium Bicarbonate Buffer (pH=8.5)
2. Transfer IgG solution into HRP,SE vial, mix well
3. Reaction for 1~2 hours at room temperature, gently shaking during the reaction is preferred.
4. Add equal volume of 100mM Tris-HCl Buffer (pH=7.5)+ 0.5% BSA
5. Store the mixture at 4°C, and it is ready to use.