

402-0005 Protocol

Maleimide Activated Alkaline Phosphatase Protocol (402-0005)

Introduction

Maleimide-activated Alkaline Phosphatase is designed to facilitate the conjugation of the Alkaline Phosphatase to proteins, peptides and ligands that contain sulfhydryl (-SH) groups. The reactive maleimide group that has been added to the Alkaline Phosphatase allows conjugation reactions to be performed very efficiently at physiological pH.

Storage of Maleimide-activated Alkaline Phosphatase

Maleimide-activated Alkaline Phosphatase is shipped at ambient temperature as a lyophilized powder plus special stabilizers and should be stored upon receipt at -20 degrees celcius. Shelf life is >1 year.

Protocol

1. Allow the vial of maleimide-activated Alkaline Phosphatase to warm to room temperature.
2. Remove the crimp seal and reconstitute the lyophilized powder by adding a buffered solution of the thiol-containing molecule. Gently agitate the vial to ensure thorough dissolution.
Note: The buffer in which the thiolated molecule is dissolved should be in the pH range 6.5-7.5. Phosphate buffered saline pH 7.2 is commonly employed.
3. Incubate for 3-4 hours (or overnight) at room temperature and then store at 4 degrees Celsius.