

408-0005 Protocol

Maleimide Activated BSA Protein Protocol (408-0005)

Introduction

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

Storage of Maleimide-activated BSA

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

Protocol

1. Allow the vial of maleimide-activated BSA 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.