

General Actin Buffer

Cat. # BSA01

Upon arrival store at 4°C (desiccated)

See datasheet for storage after reconstitution

Material

General Actin Buffer is supplied as a white lyophilized powder.

Storage and Reconstitution

Lyophilized General Actin Buffer is stable at 4°C desiccated (<10% humidity) for 1 year.

Each bottle of General Actin Buffer should be resuspended in 100 ml of sterile de-ionized water to give a 1X strength buffer; 0.2 mM calcium chloride, 5 mM Tris-HCl pH 8.0.

Uses

Used as an in vitro buffer for monomer actin proteins from a wide range of species. This is the buffer of choice when working with biologically active actin proteins. The buffer should be supplemented with 0.2 mM ATP (Cat. # BSA04-001) to create an optimal actin monomer buffer. Calcium ions and ATP are required for actin conformation, ATP is hydrolysed during actin polymerization and is required for the polymerization process.

Product Citations / Related Products

For the latest citations and related products please visit www.cytoskeleton.com.