

## General Actin Buffer

**Cat. # BSA01**

**Amount: see reconstitution**

**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 sterile de-ionized water according to the table below to give a 1X strength buffer: 0.2 mM calcium chloride, 5 mM Tris-HCl pH 8.0.

Cat #	Resuspension volume
BSA01-001	10 ml
BSA01-010	100 ml

### 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. Additionally, ATP is required for the polymerization process and is hydrolyzed during actin polymerization.

### Product Citations / Related Products

For the latest citations and related products please visit [www.cytoskeleton.com](http://www.cytoskeleton.com).