

## Product Information

### EasyProbes™ ActinGreen 488 Stain

Cat. No.	Unit Size
C060	2×2.5 mL

**Storage upon receipt:**

- 2-8°C
- Protect from light

**Ex/Em:** 500/525 nm,

### Product Description

Our actin probes are prepared by conjugating phalloidin with our Andy Fluor dyes. These fluorescently-labeled phalloidins have virtually identical binding properties with actin from different species including plants and animals. These phalloidin conjugates maintain high binding affinity and selectivity with F-actin, providing useful probes for multicolor imaging applications.

### Feature:

- **High selectivity with F-actin**
- **Multicolor selection**
- **Good photostability**
- **Superior to antibody staining**

### Stain Protocol

This procedure may not be optimum for a particular experimental system, but has yielded consistent results in most instances. The following protocol describes the staining procedure for adherent cells grown on glass coverslips.

- 1.1 Wash cells twice with prewarmed phosphate-buffered saline, pH 7.4 (PBS).
- 1.2 Fix the sample in 3.7% formaldehyde solution in PBS for 10 minutes at room temperature. **Note: Methanol can disrupt actin during the fixation process. Therefore, it is best to avoid any methanol containing fixatives. The preferred fixative is methanol-free formaldehyde.**
- 1.3 Wash two or more times with PBS.
- 1.4 Place each coverslip in a glass petri dish and extract it with a solution of acetone at  $-20^{\circ}\text{C}$  or 0.1% Triton X-100 in PBS for 3 to 5 minutes.
- 1.5 Wash two or more times with PBS.
- 1.6 Pre-incubate cells with PBS containing 1% BSA for 20–30 minutes.
- 1.7 Apply 2 drops of **EasyProbes™ ActinGreen 488 Stain** on the coverslip, and incubate for 30 minutes at room temperature. To avoid evaporation, keep the coverslips inside a covered container during the incubation.
- 1.8 Wash two or more times with PBS.
- 1.9 For long-term storage, the cells should be air dried and then mounted in a permanent mountant such as Cytoseal. Specimens prepared in this manner retain actin staining for at least six months when stored in the dark at 2–6°C.