

Instruction Manual

Product Name	Product Description	Size	Catalog Number
PromoFluor Antifade Reagent	PromoFluor Antifade Reagent	8 ml	PK-PF-AFR1

Product Description

Loss of fluorescence signals due to irreversible photobleaching processes may result in significant reduction of sensitivity, particularly when high intensity excitation light is used for a longer duration and target molecules in your sample are of low abundance.

PromoFluor Antifade Reagent minimizes photobleaching by stabilizing fluorochromes in fixed cells, fixed tissues and cell-free preparations. It is compatible with a multitude of dyes and dye complexes and very useful particularly for multicolor applications. It comes premixed and ready to use in a convenient pipet flask — just add a drop of reagent and mount. Samples can be saved for months after mounting retaining fluorescence during prolonged storage.

PromoFluor Antifade Reagent is an ideal mounting medium for immunofluorescence microscopy allowing you to get the finest images from your precious immunofluorescence preparations. It significantly increases the photostability of many common fluorophores resulting in prolonged gleaming signals and optimal contrast of the images. Moreover, it gently solidifies after the attachment of the cover slide thus permanently conserving the cell preparation and fixing the cover slide making the use of nail polish for slide immobilization superfluous.

PromoFluor Antifade Reagent is stable at 4°C in a dark place for about 4 weeks. For long-term storage it is strongly recommended to make aliquots and store the tubes tightly capped at -20°C to avoid polymerization. Before use, warm tubes to room temperature. Discard, if any crystalline material is seen in the tube or on the slide.

Protocol

- Complete staining (and counterstaining).
- If PromoFluor Antifade Reagent has been stored at -20°C make sure to thaw the reagent completely before use. It can be warmed to 37°C for a short period to thaw it properly and should be a viscous liquid then.
- Prior to mounting slides, remove excess moisture from the slide.

For mounting slides, apply one small drop of PromoFluor Antifade Reagent onto tissue sections or cell preparations which are mounted on a glass microscope slide. Then coverslip the slide and allow the PromoFluor Antifade Reagent to disperse over the entire section. (For specimens mounted on coverslips, place a drop of Antifade Reagent onto a clean slide and carefully lower the coverslip onto the Antifade Reagent). Do not apply an excess of mounting medium and avoid trapping any air bubbles.

- The coverslips should not be moved until the mounting medium has started to cure (at RT!). Evaluate slides immediately or store them at 4°C in the dark for later use.
- PromoFluor Antifade Reagent will slowly solidify at RT within a few hours (typically within 1-2 hours) starting at the edge of the coverslip. This prevents the cells from drying out and permanently immobilizes the coverslip.
- After solidification, store slides at 4°C in the dark.

To further reduce photobleaching, expose samples only when observing or recording and avoid applying too high excitation light intensity (use neutral density filters). Moreover, you may use appropriate high-quality optical filters, high-numerical aperture objectives, relatively low magnification and high-speed films or high-efficiency detectors.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.