

# VECTASHIELD®

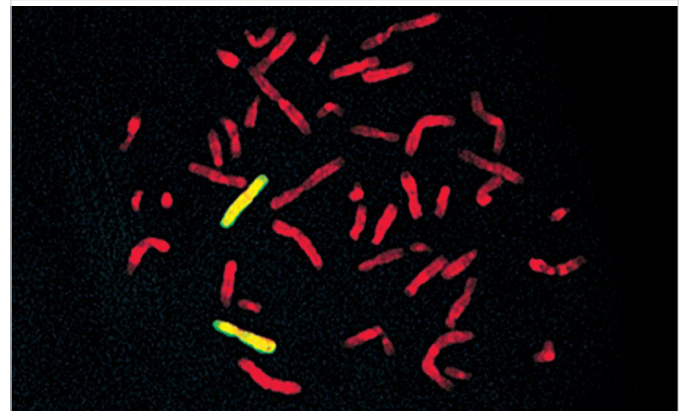
## Antifade Mounting Medium with PI

<b>Cat. No.</b>	<b>H-1300</b>
<b>Storage</b>	Refrigerate
<b>Concentration of PI</b>	1.5 µg/ml
<b>Refractive Index</b>	1.45

**Description** VECTASHIELD Antifade Mounting Medium is a unique, stable formula for preserving fluorescence. VECTASHIELD Antifade Mounting Medium prevents rapid photobleaching of fluorescent proteins and fluorescent dyes (see website for compatibility).

Each lot is tested for photobleaching retardance with qualitative and quantitative methods.

This special formulation of VECTASHIELD Antifade Mounting Medium contains propidium iodide (PI), which fluoresces when bound to DNA/RNA and is used as a nuclear or chromosomal counterstain. This product is designed to be used for in situ hybridization techniques or other procedures requiring fluorescent labeling of DNA. PI excites at about 535 nm and emits at about 615 nm when bound to DNA, producing a red fluorescence.



Human chromosome 6 paint detected with FITC-labeled probe (yellow) and mounted with VECTASHIELD Mounting Medium with PI (red). Photo courtesy of CytoCell Ltd.

### Application

To mount tissues or cells on a slide dispense VECTASHIELD Mounting Medium onto the specimen using a disposable pipet or pipet tip. Small drop volumes of approximately 25 µl (per 22 mm x 22 mm coverslip) are recommended. Coverslip the specimen and allow VECTASHIELD Mounting Medium to disperse over the entire section. Tightly seal the bottle with the screw cap after use and store in an upright position.

VECTASHIELD Antifade Mounting Medium does not solidify, but remains a liquid on the slide and can be stored without sealing. If desired, coverslips can be permanently sealed around the perimeter with nail polish or a plastic sealant. Mounted slides should be stored at 4 °C protected from light.

Note: PI is a mutagen and appropriate precautions should be taken when using it.

### VECTASHIELD Antifade Mounting Media Products:

Product	Counterstain	Cat. No.	Unit Size	Hardening
VECTASHIELD® Antifade Mounting Medium	none	H-1000	10 ml	no
	DAPI	H-1200	10 ml	no
	PI	H-1300	10 ml	no
VECTASHIELD® PLUS Antifade Mounting Medium	none	H-1900	2 ml, 10 ml	no
	DAPI	H-2000	2 ml, 10 ml	no
VECTASHIELD® HardSet™ Antifade Mounting Medium	none	H-1400	10 ml	yes
	DAPI	H-1500	10 ml	yes
	TRITC-Phalloidin	H-1600	10 ml	yes
VECTASHIELD Vibrance® Antifade Mounting Medium	none	H-1700	2 ml, 10 ml	yes
	DAPI	H-1800	2 ml, 10 ml	yes