

VECTASHIELD® HardSet™

Antifade Mounting Medium with TRITC-phalloidin

Cat. No.	H-1600
Storage	Refrigerate
Concentration of TRITC-phalloidin	0.2 µg/ml
Refractive Index	1.36 (initial) 1.46 (cured)

Description

VECTASHIELD HardSet Antifade Mounting Medium preserves fluorescence and hardens after coverslipping. VECTASHIELD HardSet Antifade Mounting Medium's unique, stable formula prevents rapid photobleaching of fluorescent proteins and fluorescent dyes (see website for compatibility).

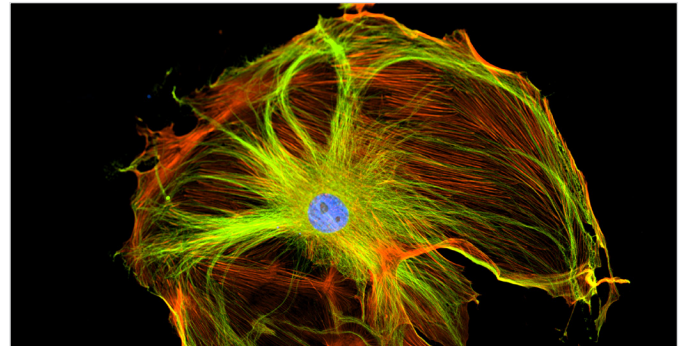
Each lot is tested for photobleaching retardance with qualitative and quantitative methods.

This special formulation of VECTASHIELD HardSet Antifade Mounting Medium contains TRITC-phalloidin. Phalloidin is a bicyclic heptapeptide that specifically binds at the interface between the F-actin subunits. Fluorescent derivatives of phalloidin are used to stain actin filaments. TRITC (tetramethylrhodamine) is excited at 544 nm and emits at 572 nm, producing an orange-red fluorescence.

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.

After approximately 15 minutes at room temperature, the coverslip will become immobilized, and optimal antifade ability and refractive index



Mouse embryonic fibroblast stained with mouse anti-tubulin followed by fluorescein horse anti-mouse IgG (green) and mounted in VECTASHIELD Hardset Antifade Mounting Medium with DAPI (blue) and TRITC-phalloidin (red).

will be achieved. After curing at room temperature for 15 minutes, slides can be placed at 4 °C, and the mounting media will harden completely overnight. Mounted slides should be stored at 4 °C or -20 °C, protected from light. For prolonged storage, -20 °C is recommended. If retraction occurs during prolonged storage, remove coverslip and remount. Coverslips can be easily removed after hardening by soaking in PBS overnight.

Notes

It is possible to get both cytoskeletal and nuclear counterstaining by mixing this mounting media with VECTASHIELD HardSet with DAPI (H-1500) at a 1:1 ratio. DAPI is a nuclear dye that is excited at 360 nm and emits at 460 nm, producing a blue fluorescence. Mix gently to avoid producing bubbles in the mixture and apply to the specimen as described above.

When mounting thick or larger sections or larger coverslips, adjustments to the mounting protocol may be required. Increase volume of mountant to provide even coverage over entire specimen extending to all edges of the coverslip.

Phalloidin is a toxin and should be handled with care.

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

For research use only. Not intended for animal or human therapeutic or diagnostic use.

©2022 Vector Laboratories. All rights reserved.

HardSet, VECTASHIELD, and VECTASHIELD Vibrance are trademarks of Vector Laboratories, Inc.
H-1600.LBL-02356.Rev00

6737 Mowry Avenue, Newark CA 94560

vectorlabs.com