

VECTASHIELD® HardSet™

Antifade Mounting Medium

Cat. No. H-1400

Storage Refrigerate

Refractive Index 1.36 (initial)
1.46 (cured)

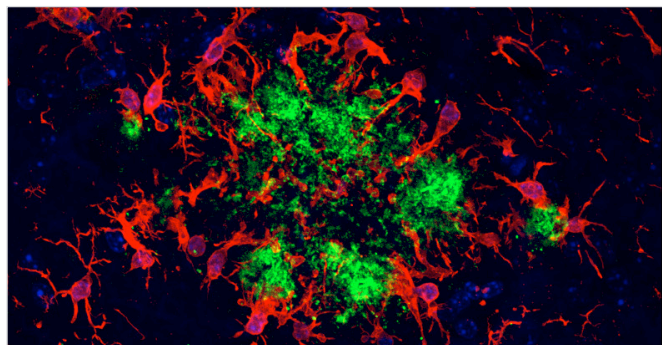
Description VECTASHIELD HardSet Antifade Mounting Medium preserves fluorescence and hardens after coverslipping. VECTASHIELD HardSet Antifade Mounting Medium's unique, stable formula prevents rapid photobleaching of fluorescent proteins and fluorescent dyes (see website for compatibility).

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

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 25µl (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 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



Alzheimer's disease mouse brain model, double label: glial cells stained with Iba1 (rp) followed by DyLight™ 594 Goat Anti-Rabbit IgG; plaques stained with Biotinylated Mouse Anti-Human Amyloid Beta followed by DyLight 488 Streptavidin. Counterstained with Hoechst 33342 and mounted with VECTASHIELD HardSet Antifade Mounting Medium.

from light. For prolonged storage, -20 °C is recommended. If retraction occurs during prolonged storage, remove coverslip and remount. Coverslips can be easily removed by soaking in PBS overnight.

Notes:

When mounting thicker or larger sections or using 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. Use of non-hardening VECTASHIELD Antifade Mounting Medium (H-1000) may be preferred.

The nuclear counterstains, DAPI and PI, will diffuse out of nuclei over time if the cells have been fixed in 100% methanol. In this case non-hardening VECTASHIELD Antifade Mounting Medium may be preferred.

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.

DyLight is a trademark of Pierce Biotechnology, Inc.

H-1400.LBL-02354.Rev00