

VECTASHIELD® PLUS

Antifade Mounting Medium with DAPI

Cat. No. H-2000

Storage 2–8 °C

Concentration of DAPI 1.0 µg/ml

Refractive Index 1.45

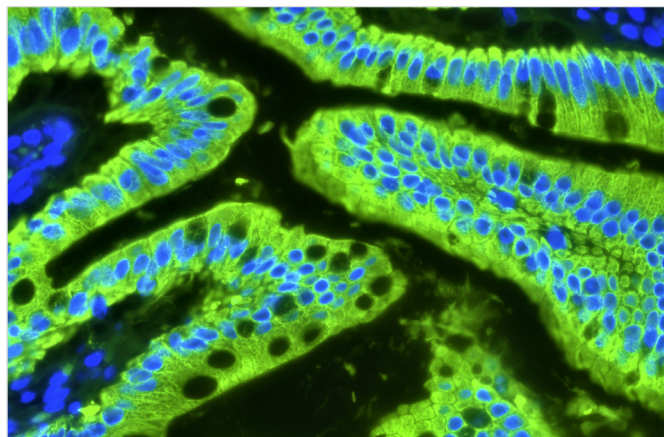
Description VECTASHIELD PLUS Antifade Mounting Medium is a unique, stable formula for preserving fluorescence. VECTASHIELD PLUS Antifade Mounting Medium prevents rapid photobleaching of fluorescent proteins and fluorescent dyes (see website for compatibility).

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

This formulation of VECTASHIELD PLUS Antifade Mounting Medium contains 4',6-diamidino-2-phenylindole (DAPI), which fluoresces when bound to DNA. DAPI excites at about 360 nm and emits at about 460 nm when bound to DNA, producing a blue fluorescence, and is ideal as a nuclear or chromosomal counterstain.

Application

To mount tissues or cells on a slide dispense VECTASHIELD PLUS Antifade Mounting Medium onto the specimen using a disposable pipet



Colon: Rabbit Anti-Cytokeratin detected with DyLight™ 488 Anti-Rabbit (green, DI-1488). Mounted in VECTASHIELD PLUS Antifade Mounting Medium with DAPI (nuclei, blue).

or pipet tip. Coverslip the specimen and allow VECTASHIELD PLUS Antifade 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 PLUS 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 2–8 °C protected from light.

Slides can be viewed 30 minutes after mounting, however optimal antifade performance is achieved after 1 hour.

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-2000.LBL-00830.Rev.00