

Thermo Scientific™ Shandon Tissue Marking Dyes Instructions for Use

Thermo Scientific Shandon Tissue Marking Dyes are intended as an aid in the orientation of excised surgical specimens through frozen sectioning, or processing and sectioning, and subsequent microscopic examination of tissues. The marking of multiple margins with these dyes is especially useful in the technique known as Mohs chemosurgery.

Instructions For Use

These dyes are normally applied to fresh tissue, but may also be used with fixed tissue. Whether fresh or fixed, the tissue should be blotted to remove excess fluid from the surface before applying the dye. The dyes may be applied using the wooden sticks provided, or, for larger surfaces, cotton swabs may be more useful.

Warning: Do not apply these dyes to a living patient. These dyes will cause permanent coloration.

Dyes will permanently bind to the tissue in 1 to 2 minutes after application. Do not apply heat. The tissue can then be frozen and prepared for sectioning, or placed into formalin or other fixatives and processed for paraffin embedding.

Order Information

Product	Size	Qty.	Ref
Tissue Marking Dyes Kit: (includes 28mL each of blue, black, yellow, green and red dyes, wooden holding block, and 50 applicator sticks.)		Ea.	3120123
Blue Tissue Marking Dye	28mL	Ea.	3120124
Black Tissue Marking Dye	28mL	Ea.	3120125
Yellow Tissue Marking Dye	28mL	Ea.	3120126
Red Tissue Marking Dye	28mL	Ea.	3120127
Green Tissue Marking Dye	28mL	Ea.	3120128

