

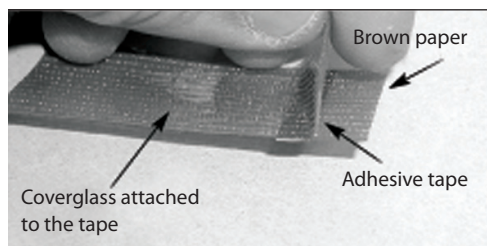
CISH UnderCover™ Slip

Cat. nos. 00-8403 and 00-8404

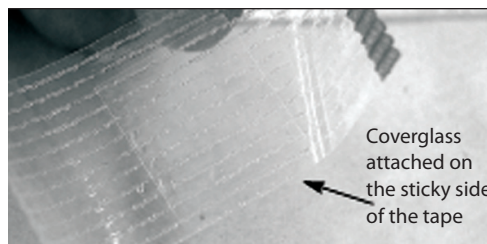
Intended use: For research use only. CAUTION: Not intended for human or animal therapeutic or diagnostic use.

Invitrogen's CISH UnderCover™ Slips consist of a coverslip attached to adhesive tape, offering a convenient alternative to traditional coverslip sealing for *in situ* hybridization reactions. Eliminate the hassle of messy rubber cement sealing under a ventilated fume hood. CISH UnderCover™ Slips form a heat-resistant, easily-applied, peel-and-stick coverslip for hybridization, sealed to prevent probe evaporation and tissue specimen dehydration on glass microscope slides. Once hybridization is complete, UnderCover™ Slips peel off cleanly and quickly. Available in 18x18 mm (Cat. no. 00-8403, Qty: 25) and 22x22 mm (Cat. no. 00-8404, Qty: 25) coverslip sizes for a variety of different sized tissue specimens.

Instructions for use:



1. Remove the tape from the brown paper backing.



2. Structure of the CISH UnderCover™ Slip.

Continued on back side.

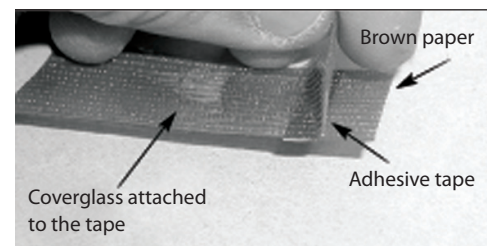
CISH UnderCover™ Slip

Cat. nos. 00-8403 and 00-8404

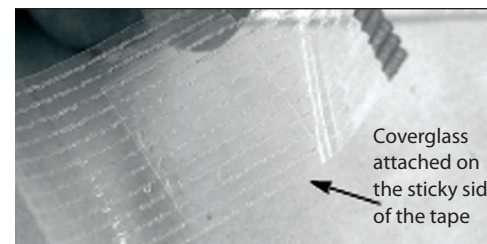
Intended use: For research use only. CAUTION: Not intended for human or animal therapeutic or diagnostic use.

Invitrogen's CISH UnderCover™ Slips consist of a coverslip attached to adhesive tape, offering a convenient alternative to traditional coverslip sealing for *in situ* hybridization reactions. Eliminate the hassle of messy rubber cement sealing under a ventilated fume hood. CISH UnderCover™ Slips form a heat-resistant, easily-applied, peel-and-stick coverslip for hybridization, sealed to prevent probe evaporation and tissue specimen dehydration on glass microscope slides. Once hybridization is complete, UnderCover™ Slips peel off cleanly and quickly. Available in 18x18 mm (Cat. no. 00-8403, Qty: 25) and 22x22 mm (Cat. no. 00-8404, Qty: 25) coverslip sizes for a variety of different sized tissue specimens.

Instructions for use:

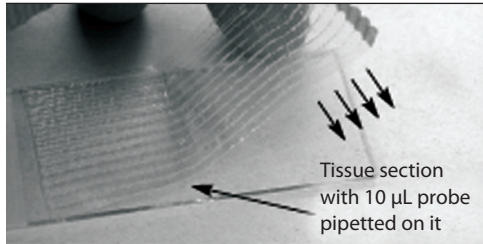


1. Remove the tape from the brown paper backing.

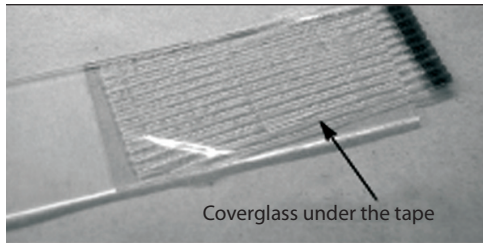


2. Structure of the CISH UnderCover™ Slip.

Continued on back side.

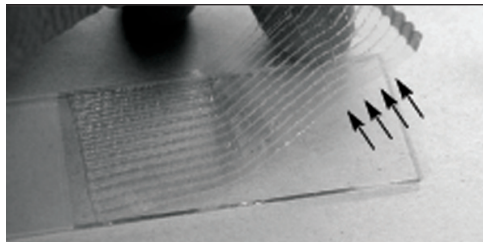


3. Apply the tape to the slide with coverglass facing the tissue section.



4. A sealed slide, ready for denaturation and hybridization.
(No humidity chamber required during hybridization)

Precaution: Do not apply extra pressure to prevent liquid from escaping under the coverglass.



5. Peel off tape after hybridization. Proceed to stringency washes and CISH or FISH detection as usual.

Storage: Store at room temperature.

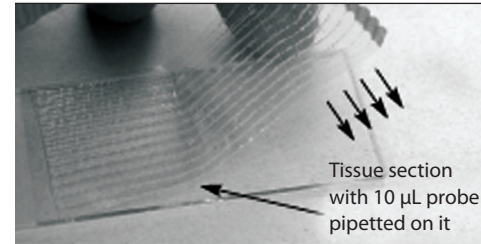
 **invitrogen**[™]

www.invitrogen.com

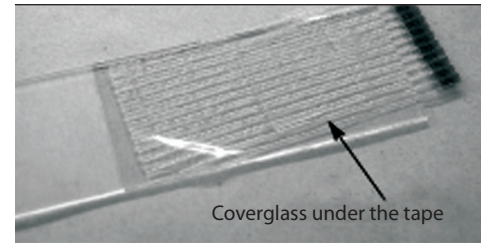
Invitrogen Corporation | 542 Flynn Road | Camarillo, CA 93012

t: 800 955 6288 | e: tech_support@invitrogen.com

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

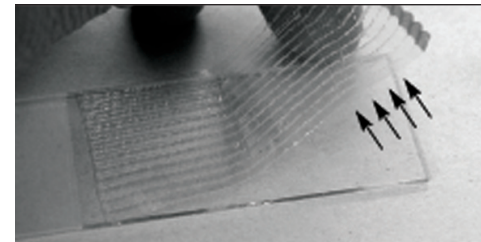


3. Apply the tape to the slide with coverglass facing the tissue section.



4. A sealed slide, ready for denaturation and hybridization.
(No humidity chamber required during hybridization)

Precaution: Do not apply extra pressure to prevent liquid from escaping under the coverglass.



5. Peel off tape after hybridization. Proceed to stringency washes and CISH or FISH detection as usual.

Storage: Store at room temperature.

 **invitrogen**[™]

www.invitrogen.com

Invitrogen Corporation | 542 Flynn Road | Camarillo, CA 93012

t: 800 955 6288 | e: tech_support@invitrogen.com

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