

IMMUNOHISTOCHEMISTRY (IHC-FR) - FROZEN SECTIONS PROTOCOL

Frozen sections: Once mounted on APES coated slides, frozen sections are best kept at -80°C until needed.

1. When required, leave to warm at room temperature for 5 min.
2. Pre-cool the fixative (acetone, methanol or ethanol) (Abcam recommends starting with acetone) at -20°C for 30 min.
3. Fix with the pre-cooled fixative for 5 -10 min, at room temperature
4. Rinse 3-4 X in PBS.
5. Continue with the immunohistochemical staining protocol. The absence of formalin eliminates the need for an antigen retrieval step. However, if frozen tissue or cytological specimens have been fixed in formalin, antigen retrieval can be attempted although the friable nature of the specimens may compromise the success. *See IHC-Paraffin protocol (IHC-P) , A) and B) for detailed protocols for chromogenic and fluorescent detection.*