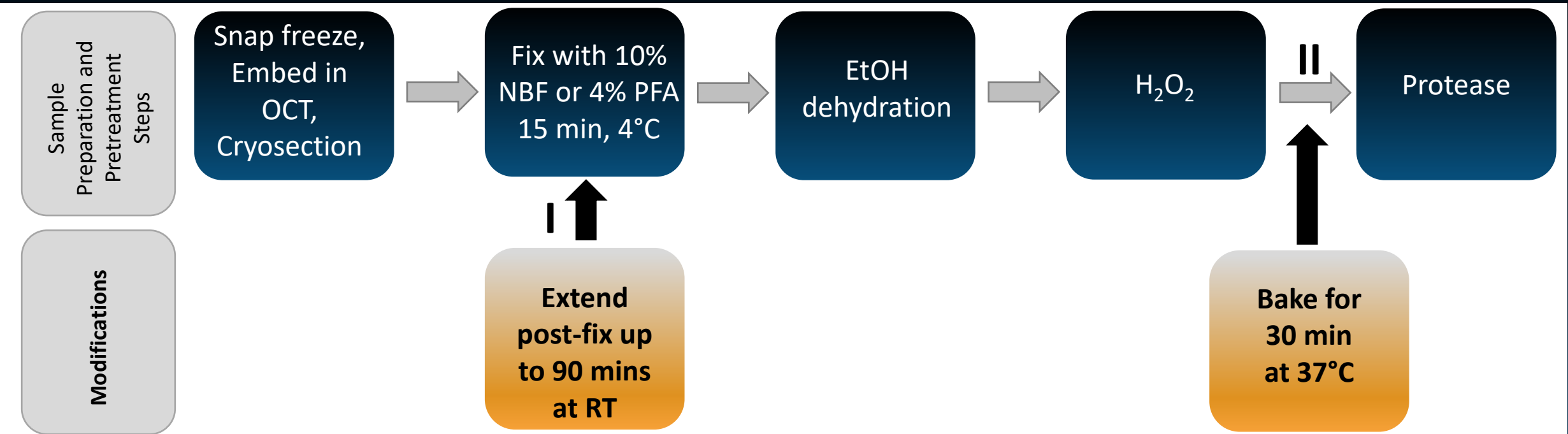


SAMPLE DETACHMENT: FRESH FROZEN SAMPLE



- Increase post-fix time and add additional baking step after H₂O₂
- Recommended section thickness 10-20 μm
- Always use Superfrost Plus slides (Fisher Scientific, Cat #12-550-15)
- Bake/Dry slides in active air circulating oven or HybEZ oven (Bake mode)