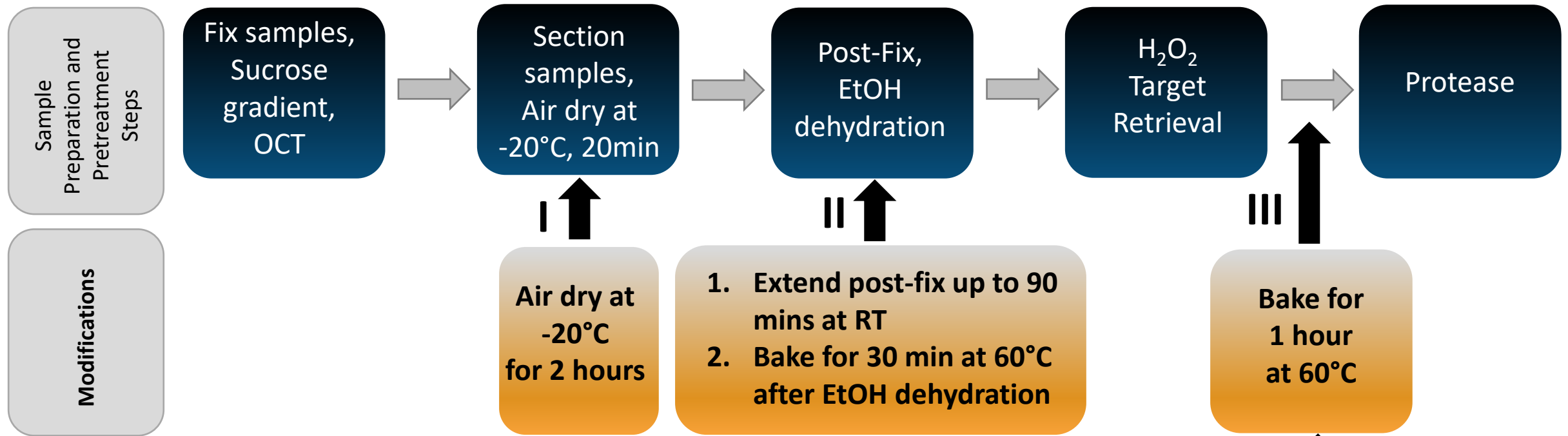


SAMPLE DETACHMENT: FIXED FROZEN SAMPLE



- Add additional steps: See I, II, and III above
- Recommended section thickness 7-15 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)
- Maintain mild boiling of target retrieval at $\sim 100^\circ\text{C}$

Omit Target Retrieval and extend protease time to 45 min if above modifications (I, II, III) fail