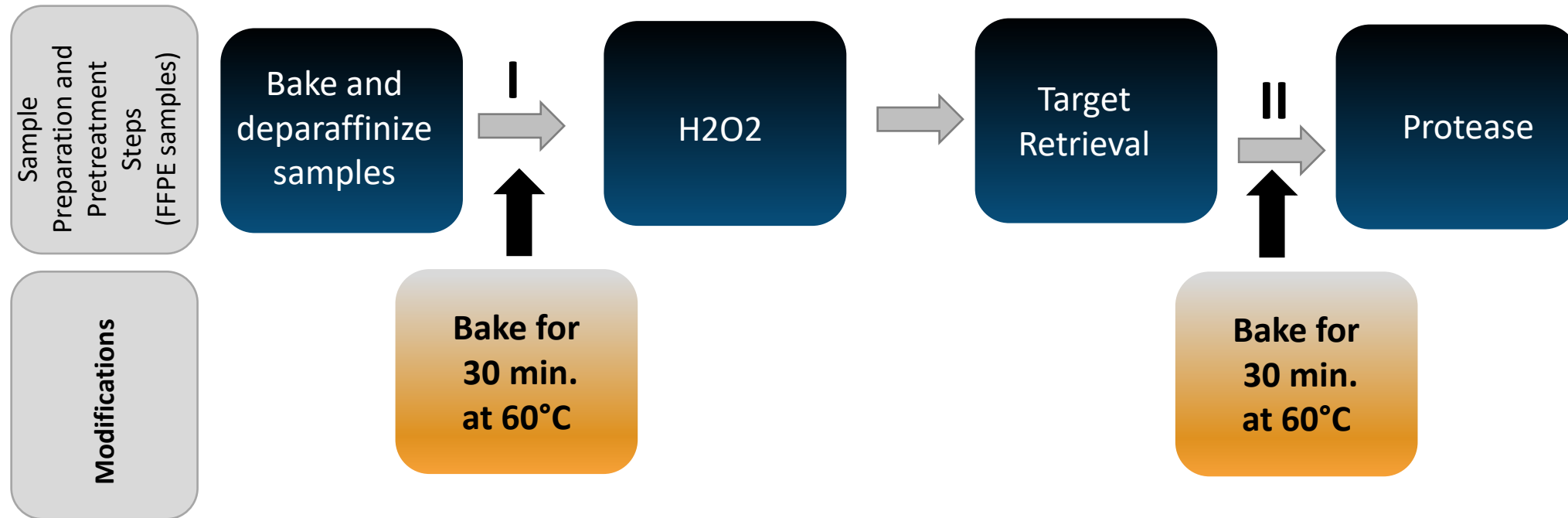


SAMPLE DETACHMENT: FFPE SAMPLES



- Add additional baking steps: See I and II above
- Always use Superfrost Plus Slides (Fisher Scientific, Cat #12-550-15)
- Bake/Dry slides in active air circulating oven
- Maintain mild boiling of target retrieval at ~100°C