



## Stripping Procedure

---

1. Put the membrane back to a hybridization tube. Add 30 ml of Stripping buffer 0.4M NaOH. Incubate with rotating at 45 °C for 30 min.
2. Decant the buffer and incubate the membrane with 30 ml 0.2M Tris-HCL, pH 7.6, 0.1X SSC, and 0.1% SDS at 45°C for 15 min.
3. Membranes are ready for hybridization or air dry the membrane at room temperature.
4. Store dry membrane protected between papers at room temperature until ready for further use.

Note: We do not encourage stripping the array membranes more than three times.