

Setting up a ligation reaction with the Quick Ligation Kit (M2200)

Overview

Protocol

1. Combine 50 ng of vector with a 3-fold molar excess of insert. Adjust volume to 10 μ l with dH₂O.
2. Add 10 μ l of 2X Quick Ligation Buffer and mix.
3. Add 1 μ l of Quick T4 DNA Ligase and mix thoroughly.
4. Centrifuge briefly and incubate at room temperature (25°C) for 5 minutes.
5. Chill on ice, then transform or store at -20°C.
6. Do not heat inactivate. Heat inactivation dramatically reduces transformation efficiency.