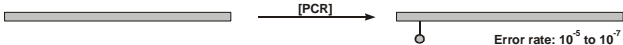


**1. Proofreading enzyme**



**2. Non-proofreading enzyme (e.g. Taq-polymerase)**



**3. Non-proofreading enzyme (e.g. Taq-polymerase)  
under error-prone conditions (e.g.  $Mn^{2+}$ )**

