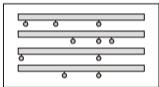


Starting pool of genes

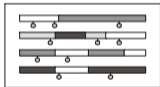
*Single gene shuffling*



*Multiple gene shuffling*



Library of point mutants



Library of chimaeras

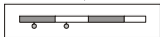
*Selection*



*Selection*



Customized gene



Customized gene