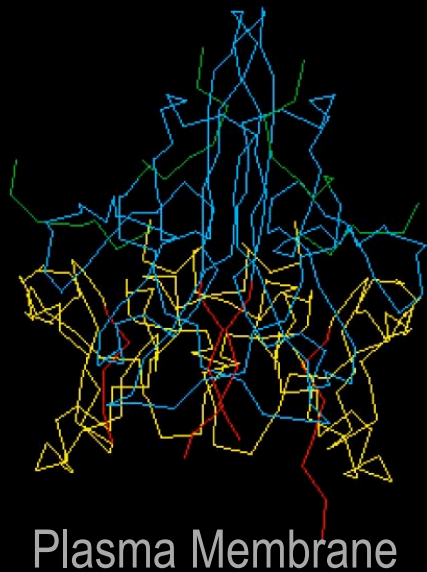
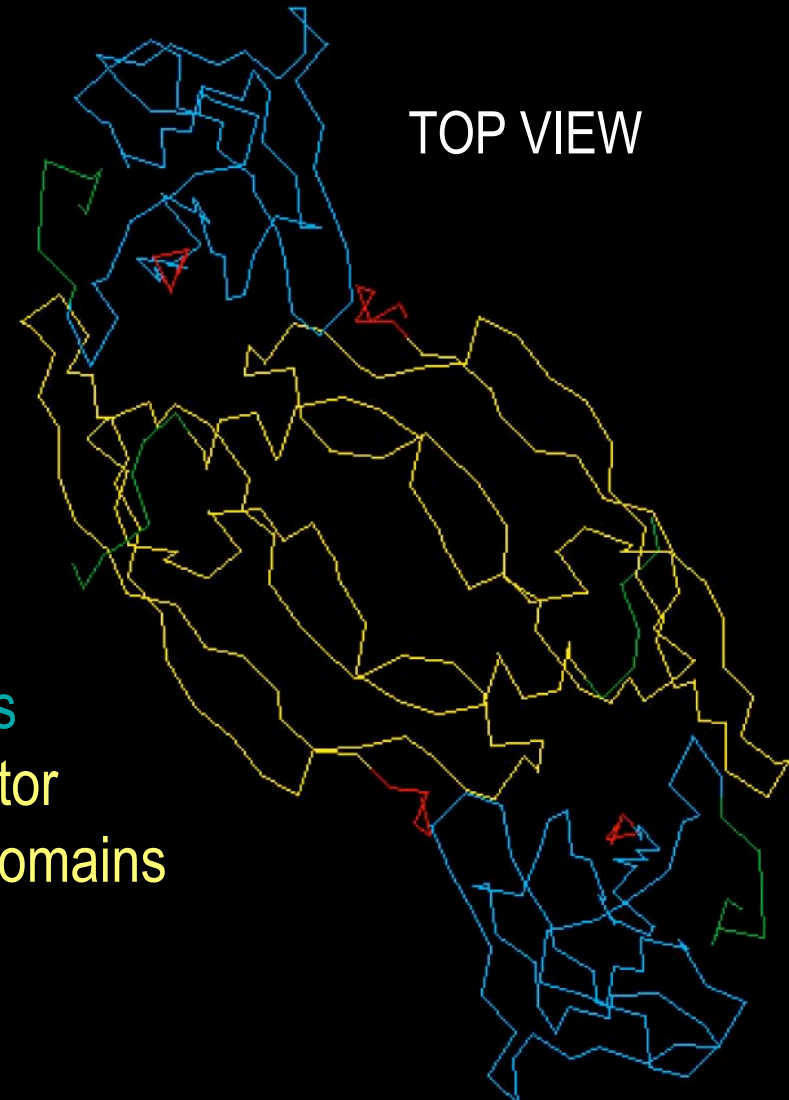
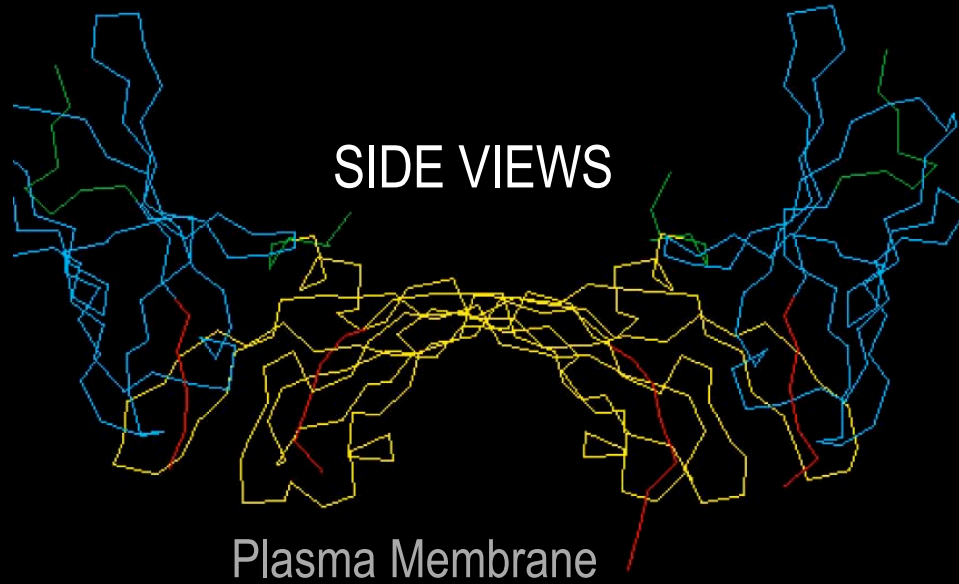


Stick Models of VEGF Binding to VEGF Receptor



- 2 VEGF chains
- 2 VEGF receptor extracellular domains
- N-termini
- C-termini