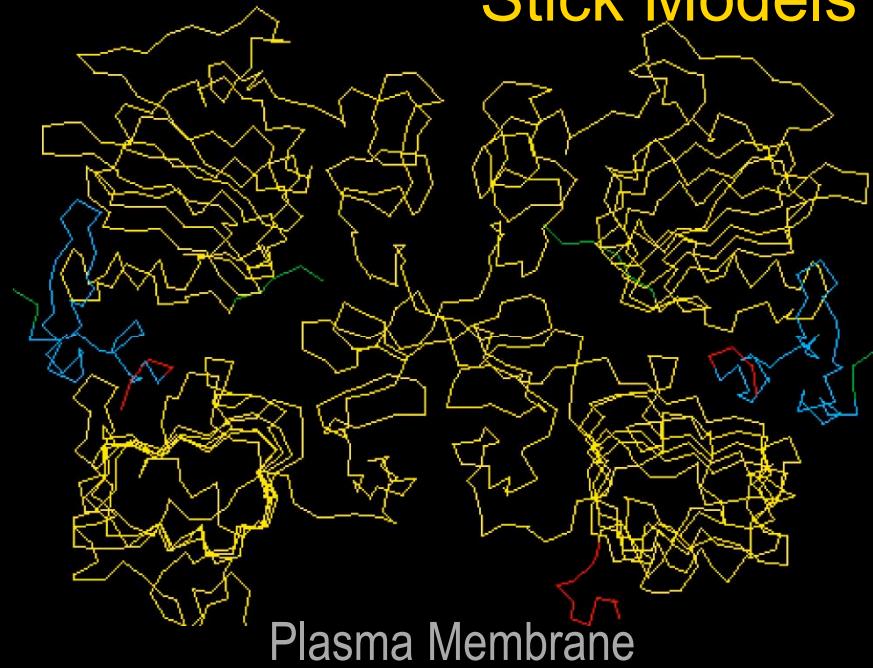


Stick Models of EGF binding to EGF Receptor Extracellular Domains



- EGF
- EGF receptor extracellular domain

SIDE VIEWS

