

Amino Acid Sequence of Proteins

- Proteins are long beaded chains of typically hundreds of amino acids.

P R C T E I N S E Q V E N C E

Working Protein (A,C,D,E,F,G,H,I,K,L,
Gene M,N,P,Q,R,S,T,V,W,Y)
Product