

SH2 and SH3 Domain Distribution in Proteins

