

Examples of Myristoylated Proteins

Src

Lck

Abl

cAMP-dependent protein kinase

Calcineurin (PP2B)

Gat (Transducin)

Gai1

Retroviral gag

Picornavirus VP4

HIV p28 Arf