

Examples of Farnesylated Proteins

Ras-H, -N, -KA + -KB

Lamins A + B

RhoA + B

Rap2

Rheb

Phosphorylase kinase $\alpha + \beta$

Rhodopsin kinase

Transducin γ

cGMP phosphodiesterase α

IFN-induced guanylate binding
protein

Ip35 phosphatase

PxF