

## Examples of Palmitoylated Proteins

Insulin receptor

HLA-DR (heavy chain)

I-A antigen

Transferrin receptor

Interleukin 1 receptor

$\beta$ -adrenergic receptor

Ankyrin

Vinculin

Nicotinic acetylcholine receptor

p21 Ras (+ other Ras  
superfamily members)

Na<sup>2+</sup> channel

G $\alpha$ s

G $\alpha$ i1

G $\alpha$ q

Myelin proteolipid protein

Rhodospin