

Examples of Glycosyl Phosphatidylinositol Anchored Proteins

Alkaline phosphatase
5' Nucleotidase
Aceylcholinesterase
Alkaline phosphodiesterase I
Variant surface glycoprotein
Thy-1
Trehalase
P63 proteinase
RT-6
Qa
ThB
T-cell activating protein
N-CAM
Heparan sulphate proteoglycan
Ornithine decarboxylase