

# 7TM Receptors Couple to G Proteins

