

Desensitization of the β -Adrenergic Receptor

