

# Light Induced Activation of cGMP Phosphodiesterase

