

# Stimuli Leading to Erk1 and Erk2 MAPK Activation

