

Conservation of Ras Signalling Pathways to Erk2 MAPK

