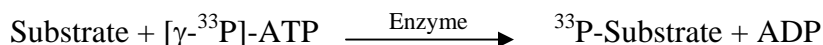


MLCK2

(Myosin light chain kinase 2-skeletal/cardiac muscle, MYLK2)

- **CAT#:** MLCK2
- **Enzyme:** Human MLCK2
- **Substrate:** Peptide substrate, [KKLNRTL^SFAEPG], 20 μM
Ca²⁺-Calmodulin, 10 μM
ATP 1 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	18.8
Ro 31-8220	660
GW 5074	2,200

- **Scientific Information:** Q9H1R3 (<http://www.expasy.org/uniprot/Q9H1R3>)

Reference compound IC50 for MLCK2/MYLK2

