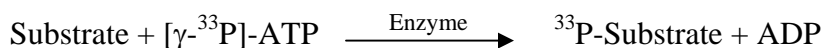


## ZAK/MLTK

(Mitogen-activated protein kinase kinase kinase MLK, MLK-like mitogen-activated protein triple kinase, Leucine zipper- and sterile alpha motif-containing kinase, Mixed lineage kinase-related kinase, Sterile alpha motif- and leucine zipper-containing kinase AZK, MLK-related kinase, MRK, Cervical cancer suppressor gene 4 protein, HCCS-4)

- **CAT#:** ZAK/MLTK
- **Enzyme:** Human ZAK/MLTK
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	38,000
Ro 31-8220	36,000
GW5074	3,500

- **Scientific Information:** Q9NYL2 (<http://www.expasy.org/uniprot/Q9NYL2>)

### Reference compound IC50 for ZAK/MLTK

