

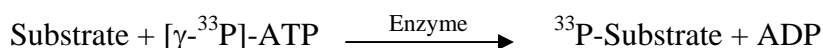
MKK6

(Constitutively active mutant S207E/T211E)

(Dual specificity mitogen-activated protein kinase kinase 6, MAP kinase kinase 6, MAPKK 6, MAPK/ERK kinase 6, SAPKK3, MAP2K6, MEK6, PRKMK6)

- **CAT#:** MKK6
- **Enzyme:** Human MKK6
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	80.0
Ro 31-8220	22,300
GW5074	46,000

- **Scientific Information:** P52564 (<http://www.expasy.org/uniprot/P52564>)

Reference compound IC50 for MKK6

