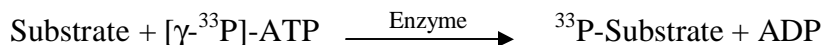


MNK2

(MAP kinase-interacting serine/threonine-protein kinase 2,
MAP kinase signal-integrating kinase 2, Mnk2, MKNK2)

- **CAT#:** MNK2
- **Enzyme:** Human MNK2
- **Substrate:** MBP (Myelin Basic Protein), 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 4.7 |
| Ro 31-8220 | 3,900 |
| GW 5074 | 1,100 |
| H-89 | 14,200 |
| Rottlerin | 556 |

- **Scientific Information:** Q9HBH9 (<http://www.expasy.org/uniprot/Q9HBH9>)

Reference Compound IC50 for MNK2

