

## P38 $\alpha$ (T106M)

(Mitogen-activated protein kinase 14, Mitogen-activated protein kinase p38 alpha, MAP kinase p38 alpha, MAP kinase MXI2, MK14)

- **CAT#:** P38 $\alpha$  (T106M)
- **Enzyme:** Human P38 $\alpha$  (T106M)
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 8,400     |
| Ro 31-8220    | 25,800    |
| GW 5074       | 18,000    |

- **Scientific Information:** Q16539 (<http://www.expasy.org/uniprot/Q16539>)

### Reference Compound IC50 for P38 $\alpha$ (T106M)

