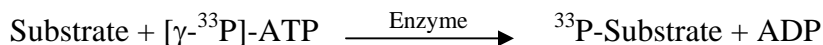


IRAK4

(Interleukin-1 receptor-associated kinase 4, IRAK-4)

- **CAT#:** IRAK4
- **Enzyme:** Human IRAK4
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.6
Ro 31-8220	950
GW 5074	10,000
H-89	>20,000

- **Scientific Information:** Q9NWZ3 (<http://www.expasy.org/uniprot/Q9NWZ3>)

Reference compound IC50 for IRAK4

