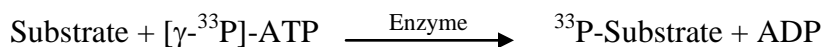


ULK2

(Serine/threonine-protein kinase ULK2, Unc-51-like kinase 2, KIAA0623)

- **CAT#:** ULK2
- **Enzyme:** Human ULK2
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	20.5
Ro 31-8220	28,600

- **Scientific Information:** Q8IYT8 (<http://www.expasy.org/uniprot/Q8IYT8>)

Reference compound IC50 for ULK2

