

MARK4

(MAP/microtubule affinity-regulating kinase 4,
MAP/microtubule affinity-regulating kinase-like 1, KIAA1860, MARKL1)

- **CAT#:** MARK4
- **Enzyme:** Human MARK4
- **Substrate:** Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 20 μ M
ATP 10 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.5
Ro 31-8220	3,300
GW5074	5,000
H-89	15,000

- **Scientific Information:** Q96L34 (<http://www.expasy.org/uniprot/Q96L34>)

Reference Compound IC50 for MARK4

