

MARK2/PAR-1Ba

(Serine/threonine-protein kinase MARK2, MAP/microtubule affinity-regulating kinase 2, ELKL motif kinase, EMK1, PAR1 homolog)

- **CAT#:** MARK2/PAR-1Ba
- **Enzyme:** Human MARK2/PAR-1Ba
- **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	1,800
GW5074	3,300
H-89	2,300

- **Scientific Information:** Q7KZ17-4 (<http://www.expasy.org/uniprot/Q7KZ17-4>)

Reference compound IC50 for MARK2/PAR-1Ba

