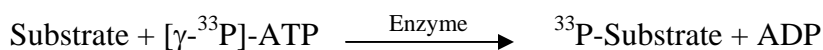


## MRCKb/CDC42BPB

(Serine/threonine-protein kinase MRCK beta, Myotonic dystrophy protein kinase-like beta, Myotonic dystrophy kinase-related CDC42-binding kinase beta, MRCK beta, DMPK-like beta, CDC42BPB)

- **CAT#:** MRCKb/CDC42BP
- **Enzyme:** Human MRCKb/CDC42BP
- **Substrate:** Peptide substrate, [KEAKEKRQEIQAKRRRLSSLRASTSKSGGSQK], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	9.8
Ro 31-8220	1,100

- **Scientific Information:** Q9Y5S2 (<http://www.expasy.org/uniprot/Q9Y5S2>)

### Reference Compound IC50 for MRCKb/CDC42BPB

