

MAPKAP kinase 3

(MAP kinase-activated protein kinase 3, MAPK3, MAPKAPK-3,
Chromosome 3p kinase, 3pK)

- **CAT#:** MAPKAPK3
- **Enzyme:** Human MAPKAPK3
- **Substrate:** Peptide substrate, [PLSRTLSTVAACK-NH₂], 5 μM
ATP 1 μM
- **Reaction:**
$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC₅₀s:**

Compounds	IC ₅₀ (nM)
Staurosporine	6,400
Ro 31-8220	31,300
GW 5074	16,800
H-89	27,200
Rottlerin	64,000

- **Scientific Information:** Q16644 (<http://www.expasy.org/uniprot/Q16644>)

Reference compound IC₅₀ for MAPKAPKinase3

