

DCAMKL1

(Serine/threonine-protein kinase DCLK1, Doublecortin domain-containing protein 3A, Doublecortin-like and CAM kinase-like 1, Doublecortin-like kinase 1, DCLK1, DCDC3A, KIAA0369)

- **CAT#:** DCAMKL1
- **Enzyme:** Human DCAMKL 1
- **Substrate:** Peptide substrate, [KKALRRQETVDAL], 20 μ M
Ca²⁺-Calmodulin, 1 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	1,750
Ro 31-8220	>20,000
GW5074	15,000
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** O15075 (<http://www.expasy.org/uniprot/O15075>)

Reference compound IC50 for DCAMKL1

