

CDK5/p35

(Cell division protein kinase 5, Tau protein kinase II catalytic subunit, TPKII catalytic subunit, Serine/threonine-protein kinase PSSALRE
P35: Cyclin-dependent kinase 5 activator 1 [Precursor], CDK5 activator 1, Cyclin-dependent kinase 5 regulatory subunit 1, Tau protein kinase II 23 kDa subunit, p23, TPKII regulatory subunit, Cyclin-dependent kinase 5 activator 1, CDK5R1)

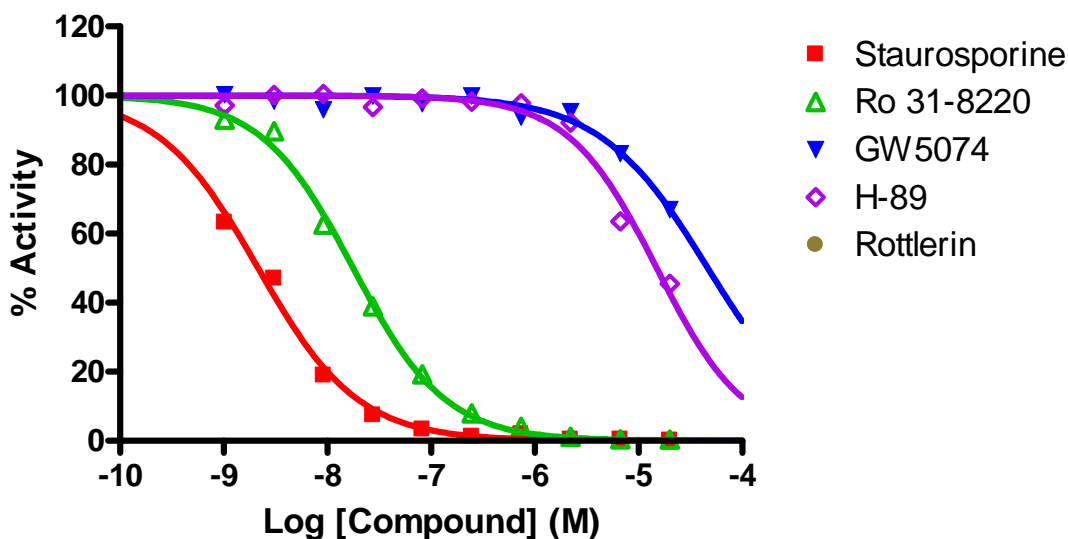
- **CAT#:** CDK5/p35
- **Enzyme:** Human CDK5/p35
- **Substrate:** Histone H1, 20 μ M
ATP 10 μ M
- **Reaction:**

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

| Compounds | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 2.1 |
| Ro 31-8220 | 18 |
| GW 5074 | >20,000 |
| H-89 | 15,000 |

- **Scientific Information:** Q00535 (<http://www.expasy.org/uniprot/Q00535>)

Reference compound IC50 for CDK5/p35



Service performed using kinases under license from  invitrogen™ part of life technologies™