

EEF2K

(Elongation factor 2 kinase, eEF-2 kinase,
Calcium/calmodulin-dependent eukaryotic elongation factor 2 kinase)

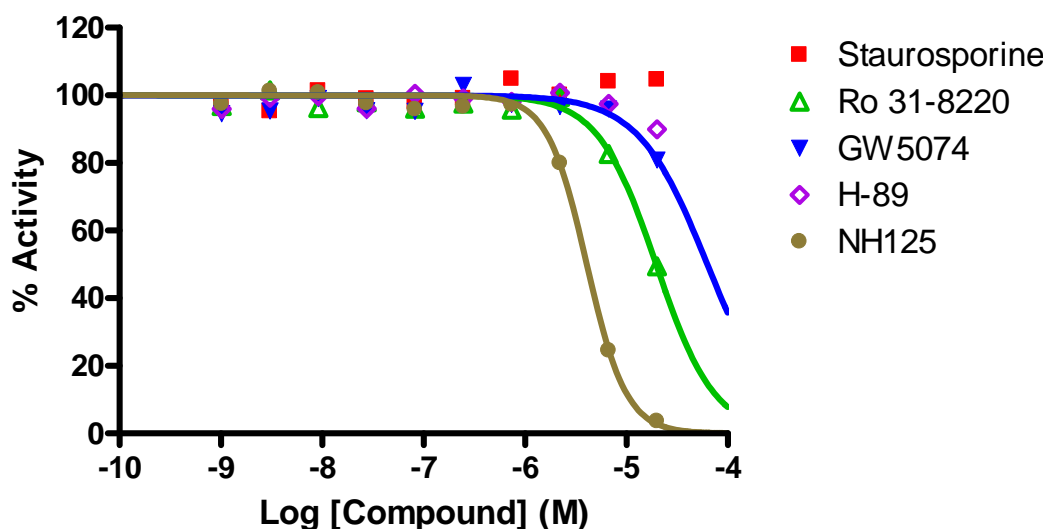
- **CAT#:** EEF2K
- **Enzyme:** Human EEF2K
- **Substrate:** Peptide substrate, [RKKFGESEKTKTKEFL], 20 μ M
Ca²⁺-Calmodulin, 1 μ M
ATP 10 μ M
- **Reaction:**

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

Compounds	IC50 (nM)
Staurosporine	>100,000
Ro 31-8220	20,000
GW 5074	63,000
NH125	4,100

- **Scientific Information:** O00418 (<http://www.expasy.org/uniprot/O00418>)

Reference compound IC50 for EEF2K



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