

## SNARK

(NUAK family SNF1-like kinase 2, SNF1/AMP kinase-related kinase, NUA2)

- **CAT#:** SNARK
- **Enzyme:** Human SNARK
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.4
Ro 31-8220	50.3
GW5074	9,900
H-89	2,000
Rottlerin	46,800

- **Scientific Information:** Q9H093 (<http://www.expasy.org/uniprot/Q9H093>)

### Reference compound IC50 for SNARK/NUAK2

