

STK16

(Serine/threonine-protein kinase 16, Protein kinase PKL12, Myristoylated and palmitoylated serine/threonine-protein kinase, MPSK, TGF-beta-stimulated factor 1, TSF-1, hPSK, MPSK1, PKL12, TSF1)

- **CAT#:** STK16
- **Enzyme:** Human STK16
- **Substrate:** MBP (Myelin Basic Protein), 10 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	770
Ro 31-8220	7,000
GW 5074	42,000
Rottlerin	175,000

- **Scientific Information:** O75716 (<http://www.expasy.org/uniprot/O75716>)

Reference Compound IC50 for STK16 10uM

