

LKB1

(LKB1: Serine/threonine-protein kinase 11, Serine/threonine-protein kinase LKB1
Renal carcinoma antigen NY-REN-19, STK11, PJS'

STRAD α : STE20-related kinase adapter protein alpha, LYK5, STRAD

MO25 α : Calcium-binding protein 39, CAB39)

- **CAT#:** LKB1
- **Enzyme:** Human LKB1 complex with STRAD α / MO25 α
- **Substrate:** Peptide substrate, [LSNLYHQGKFLQTFCSPLYRRR], 20 μ M
ATP 1 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	35.1
Ro 31-8220	3,200
GW5074	10,300
H-89	8,000

- **Scientific Information:** LKB1: Q15831 (<http://www.expasy.org/uniprot/Q15831>)
STRADA:Q7RTN6 (<http://www.expasy.org/uniprot/Q7RTN6>)
CAB39:Q9Y376 (<http://www.expasy.org/uniprot/Q9Y376>)

Reference Compound IC50 for LKB1

