
MLK3(MAP3K11)

Product code: 09-017

Human MLK3, catalytic domain [99-398 amino acids of accession number NP_002410.1] was expressed as N-terminal His-tagged protein (37kDa) using baculovirus expression system.

| | |
|-----------------------|-----------------------------|
| Assay platform | : MSA |
| Substrate | : MAP2K1 Cascade Assay * |
| ATP (μ M) | : 1000 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at 1 mM ATP (nM) | : 4.8 |

* MAP2K1/Erk2/Modified Erktide
