
TNK1

Product code: 08-104

Human TNK1, catalytic domain [106-390 amino acids of accession number NP_003976.2] was expressed as N-terminal GST-fusion protein (58 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : CSKtide |
| ATP (μ M) Km app / Bin | : 71 / 75 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 0.55 |
| IC50 at 1 mM ATP (nM) | : 1.7 |

