
HIPK1

Product code: 04-135

Human HIPK1, catalytic domain [158-555 amino acids of accession number NP_689909.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : DYRKtide-F |
| ATP (μ M) Km app / Bin | : 4.4 / 5 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 570 |
| IC50 at 1 mM ATP (nM) | : n.a. |

