
SGK3(SGKL)

Product code: 01-160

Truncated human SGK3 [119-496(end) amino acids and S486D of accession number NP_037389.4] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : SGKtide |
| ATP (μ M) Km app / Bin | : 17 / 25 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 42 |
| IC50 at 1 mM ATP (nM) | : n.a. |