

---

# AKT1

Product code: 01-101

Human AKT1, catalytic domain [104-480(end) amino acids and G478S of accession number NP\_005154.2] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP\_002604.1] using baculovirus expression system. GST-AKT1 was purified by using glutathione sepharose chromatography.

|                             |                 |
|-----------------------------|-----------------|
| Assay platform              | : MSA           |
| Substrate                   | : Crosstide     |
| ATP ( $\mu$ M) Km app / Bin | : 31 / 50       |
| Metal                       | : Mg            |
| Reference compound          | : Staurosporine |
| IC50 at ATP Bin (nM)        | : 2.7           |
| IC50 at 1 mM ATP (nM)       | : 22            |

