
PKC α (PRKCA)

Product code: 01-133

Full-length human PKC α [1-672(end) amino acids and V568I of accession number NP_002728.2] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : PKC peptide |
| ATP (μ M) Km app / Bin | : 36 / 50 |
| Metal | : Mg+Ca |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 0.33 |
| IC50 at 1 mM ATP (nM) | : 3.6 |

