
PKC γ (PRKCG)

Product code: 01-137

Full-length human PKC γ [1-697(end) amino acids of accession number NP_002730.1] was expressed as N-terminal GST-fusion protein (106 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : PKC peptide |
| ATP (μ M) Km app / Bin | : 74 / 75 |
| Metal | : Mg+Ca |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 1.1 |
| IC50 at 1 mM ATP (nM) | : 11 |

