

---

---

## PKC $\beta$ 2(PRKCB2)

Product code: 01-165

Full-length human PKC $\beta$ 2 [1-673(end) amino acids of accession number NP\_002729.2] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system.

|                             |                 |
|-----------------------------|-----------------|
| Assay platform              | : MSA           |
| Substrate                   | : PKC peptide   |
| ATP ( $\mu$ M) Km app / Bin | : 41 / 50       |
| Metal                       | : Mg+Ca         |
| Reference compound          | : Staurosporine |
| IC50 at ATP Bin (nM)        | : 0.43          |
| IC50 at 1 mM ATP (nM)       | : n.a.          |

