
CK1 γ 3(CSNK1G3)

Product code: 03-107

Full-length human CK1 γ 3 [1-447(end) amino acids of accession number NP_004375.2] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system.

| | |
|-----------------------------|--------------------|
| Assay platform | : MSA |
| Substrate | : CKtide |
| ATP (μ M) Km app / Bin | : 3.2 / 5 |
| Metal | : Mg |
| Reference compound | : 5-Iodotubercidin |
| IC50 at ATP Bin (nM) | : 920 |
| IC50 at 1 mM ATP (nM) | : >10000 |

