
CGK2(PRKG2)

Product code: 01-143

Full-length human CGK2 [1-762(end) amino acids of accession number NP_006250.1] was expressed as N-terminal GST-fusion protein (114 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.88
IC50 at 1 mM ATP (nM)	: n.a.

