
FRK

Product code: 08-167

Human FRK, catalytic domain [223-505(end) amino acids of accession number NP_002022.1] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 62 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.4
IC50 at 1 mM ATP (nM)	: 40

