
YES(YES1)[T348I]

Product code 08-533

Full-length human YES [1-543(end) amino acids and T348I of accession number NP_005424.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-YES[T348I] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 8.5 / 10
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 1.4
IC₅₀ at 1 mM ATP (nM) : 45
