
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.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 8.5 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.4
IC50 at 1 mM ATP (nM)	: 45