
TYRO3

Product code 08-109

Human TYRO3, cytoplasmic domain of [453-890(end) amino acids of accession number NP_006284.2] was expressed as N-terminal GST fusion protein (76 kDa) using baculovirus expression system. GST-TYRO3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : CSKtide

ATP (μ M) Km app / Bin : 80 / 75
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.3
IC50 at 1 mM ATP (nM) : 2.9
