
HER4(ERBB4)

Product code 08-118

Human HER4, cytoplasmic domain [676-1308(end) amino acids of accession number NP_005226.1] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system. GST-HER4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Src tide

ATP (μ M) K_{mapp} / B_{in} : 27 / 25
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 34
IC₅₀ at 1 mM ATP (nM) : 1000
