
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.

Assay platform : MSA
Substrate : Srctide

ATP (μ M) Km app / Bin : 27 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 34
IC50 at 1 mM ATP (nM) : 1000
