
HER2(ERBB2)

Product code: 08-016

Human HER2, cytoplasmic domain [676-1255(end) amino acids and P1170A of accession number NP_004439.2] was expressed as N-terminal His-tagged protein (67 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Srctide
ATP (μ M) Km app / Bin : 3.5 / 5
Metal : Mn
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 90
IC50 at 1 mM ATP (nM) : 6500
