
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) | : >10000 |

