

---

---

## 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 ( $\mu$ 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

