

## Product Information

### HER2(ERBB2)

 Product Number : **08-016**

#### Product description

Human HER2, cytoplasmic domain [676-1255(end) amino acids of accession number NP\_004439.1] was expressed as N-terminal His-tagged protein (67 kDa) using baculovirus expression system. His-tagged HER2 was purified by using Ni-NTA affinity chromatography.

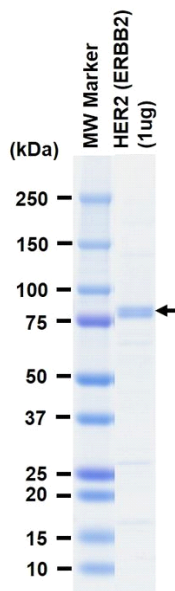
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
 5 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

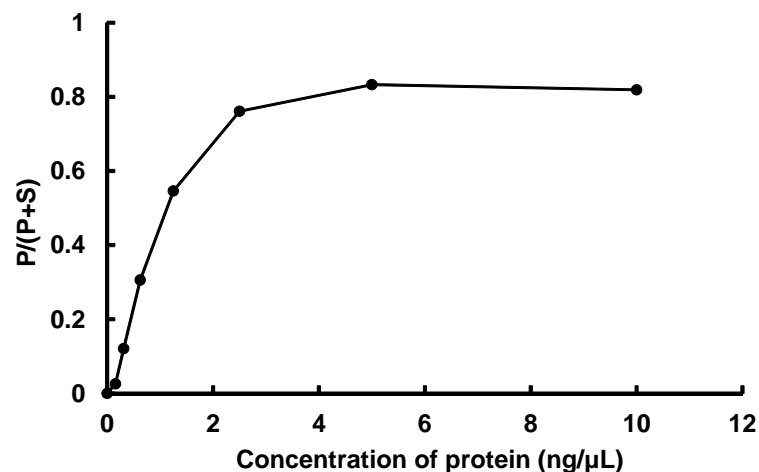
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 70 %  
 The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mn/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Srctide

ATP : 100 μM