

---

## ***EGFR***

Product code 08-115

Human EGFR, cytoplasmic domain [669-1210(end) amino acids of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP ( $\mu\text{M}$ )  $K_{\text{mapp}}$  /  $B_{\text{in}}$  : 2.7 / 5

Metal : Mg+Mn

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{\text{in}}$  (nM) : 53

IC<sub>50</sub> at 1 mM ATP (nM) : 7700

---