

---

# 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.

Assay platform : MSA  
Substrate : Srctide  
  
ATP ( $\mu$ M) Km app / Bin : 2.7 / 5  
Metal : Mg+Mn  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 53  
IC50 at 1 mM ATP (nM) : 7700

---