

---

## EGFR[d746-750/C797S]

Product code 08-564

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and C797S of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-EGFR[d746-750aa/C797S] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 8.2 / 10

Metal : Mg+Mn

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 8.0

IC<sub>50</sub> at 1 mM ATP (nM) : 130

---