

---

## EGFR[C797S/L858R]

Product code: 08-563

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and C797S/L858R of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (89kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Srctide

ATP ( $\mu$ M) Km app / Bin : 4.1 / 5

Metal : Mg+Mn

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 8.8

IC50 at 1 mM ATP (nM) : 270

---