
EGFR [T790M/L858R]

Product code 08-510

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and T790M/L858R of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[T790M/L858R] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP (μ M) K_{mapp} / B_{in} : 1.9 / 2

Metal : Mg+Mn

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 1.0

IC₅₀ at 1 mM ATP (nM) : 56
