
EGFR[d746-750]

Product code 08-527

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-EGFR[d746-750aa] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 19 / 25

Metal : Mg+Mn

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 13

IC₅₀ at 1 mM ATP (nM) : 93
