
EGFR[d746-750/T790M]

Product code: 08-528

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and T790M of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 5.4 / 5

Metal : Mg+Mn

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.52

IC₅₀ at 1 mM ATP (nM) : 9.7
