
EGFR[L861Q]

Product code: 08-513

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and L861Q of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Srctide

ATP (μ M) Km app / Bin : 7.5 / 10
Metal : Mg+Mn
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 68
IC50 at 1 mM ATP (nM) : 2200
