
EGFR[L858R]

Product code: 08-502

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and L858R 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 : 9.8 / 10

Metal : Mg+Mn

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 11

IC50 at 1 mM ATP (nM) : 360
