
EGFR

Product code: 08-115

Human EGFR, cytoplasmic domain [669-1210(end) amino acids 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 : 2.7 / 5

Metal : Mg+Mn

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 53

IC50 at 1 mM ATP (nM) : 7700
