
EGFR[D770_N771insNPG]

Product code: 08-553

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and D770_N771insNPG 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	: 2.3 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 18
IC50 at 1 mM ATP (nM)	: 930