
EGFR[T790M/L858R]

Product code: 08-510

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and T790M/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	: 1.9 / 2
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 56

