
EGFR[d746-750/T790M/C797S]

Product code: 08-565

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and T790M/C797S of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 1.8 / 2
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.84
IC50 at 1 mM ATP (nM)	: 14

