
EGFR[L861Q]

Product code: 08-513

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and L861Q 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 | : 7.5 / 10 |
| Metal | : Mg+Mn |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 68 |
| IC50 at 1 mM ATP (nM) | : 2200 |

