
FGFR2[V564I]

Product code: 08-546

Human FGFR2, cytoplasmic domain [399-821(end) amino acids and V564I of accession number NP_000132.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : CSKtide |
| ATP (μ M) Km app / Bin | : 21 / 25 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 2.0 |
| IC50 at 1 mM ATP (nM) | : 47 |
