
EPHB2

Product code: 08-129

Human EPHB2, cytoplasmic domain [581-987(end) amino acids of accession number NP_004433.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

| | |
|-----------------------------|-----------------|
| Assay platform | : MSA |
| Substrate | : Blk/Lyntide |
| ATP (μ M) Km app / Bin | : 86 / 100 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 73 |
| IC50 at 1 mM ATP (nM) | : 400 |

