
ALK[G1202R]

Product code: 08-544

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and G1202R, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 31 / 50

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 7.3

IC₅₀ at 1 mM ATP (nM) : 69
