
ALK[G1269A]

Product code 08-537

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and G1269A, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[G1269A] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Srctide

ATP (μ M) Km app / Bin : 27 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.36

IC50 at 1 mM ATP (nM) : 1.6
