
ALK[R1275Q]

Product code 08-520

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and R1275Q, 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[R1275Q] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 84 / 100

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 3.3

IC₅₀ at 1 mM ATP (nM) : 16
