
ALK

Product code 08-518

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

Assay platform : Mobility Shift Assay

Substrate : Srctide

ATP (μ M) Km app / Bin : 57 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.5

IC50 at 1 mM ATP (nM) : 15
