
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.

Assay platform : MSA
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
