
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.

Assay platform : MSA
Substrate : Srctide

ATP (μ M) Km app / Bin : 84 / 100
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 3.3
IC50 at 1 mM ATP (nM) : 16
