

---

# 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.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ 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

