

---

# ALK[F1174L]

Product code: 08-519

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and F1174L, 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	: Src tide
ATP ( $\mu$ M) $K_m$ app / $B_{in}$	: 49 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC <sub>50</sub> at ATP $B_{in}$ (nM)	: 2.4
IC <sub>50</sub> at 1 mM ATP (nM)	: 21