

---

---

# ALK[T1151\_L1152insT]

Product code: 08-539

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and T1151\_L1152insT, 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 : 110 / 100  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 6.5  
IC50 at 1 mM ATP (nM) : 16

---

---