

---

# NPM1-ALK

Product code: 08-517

Fused gene of human fusion NPM1-ALK [1-680 amino acids of accession number BAA08343.1] was expressed as N-terminal GST-fusion protein (103kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 57 / 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) : 14

---