

---

# ITK

Product code: 08-181

Full-length human ITK [2-620(end) amino acids of accession number NP\_005537.3] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 6.1 / 10

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 3.2

IC<sub>50</sub> at 1 mM ATP (nM) : 200

---