

---

## ***EPHA3***

Product code 08-122

Human EPHA3, cytoplasmic domain [579-983(end) amino acids of accession number NP\_005224.2] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-EPHA3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 170 / 150

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 37

IC<sub>50</sub> at 1 mM ATP (nM) : 76

---