

---

---

## FAK(PTK2)

Product code: 08-137

Truncated human FAK[376-1052(end) amino acids of accession number NP\_722560.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP ( $\mu$ M) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 47
IC50 at 1 mM ATP (nM)	: 230

