

---

---

## 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           |

