
IRAK1

Product code: 09-101

Truncated human IRAK1 [194-712(end) amino acids of accession number NP_001560.2] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system.

| | |
|--------------------------------|-----------------|
| Assay platform | : IMAP |
| Substrate | : SRPKtide |
| ATP (μ M) K_m app / Bin | : 27 / 25 |
| Metal | : Mg |
| Reference compound | : Staurosporine |
| IC50 at ATP Bin (nM) | : 54 |
| IC50 at 1 mM ATP (nM) | : n.a. |

