
FER

Product code: 08-139

Full-length human FER [1-822(end) amino acids of accession number NP_005237.2] was expressed as N-terminal GST-fusion protein (122 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 26 / 25

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 1.0

IC₅₀ at 1 mM ATP (nM) : 12
