
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
IC ₅₀ at ATP B_{in} (nM)	: 1.0
IC ₅₀ at 1 mM ATP (nM)	: 12