
FER

Product code: 08-139

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

Assay platform	: MSA
Substrate	: Srctide
ATP (μ M) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 12

