
CRIK(CIT)

Product code: 01-104

Human citron kinase (CRIK), catalytic domain [1-449 amino acids of accession number NP_009105.1] was expressed as N-terminal GST fusion protein (77 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Histone H3 peptide
ATP (μ M) Km app / Bin	: 7.8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 31
IC50 at 1 mM ATP (nM)	: n.a.
