
EEF2K

Product code: 10-113

Full-length human EEF2K [1-725(end) amino acids and Q361R of accession number NP_037434.2] was expressed as N-terminal GST-fusion protein (109 kDa) using E. coli expression system.

Assay platform	: MSA
Substrate	: EEF2Ktide
ATP (μ M) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: A-484954
IC50 at ATP Bin (nM)	: 330
IC50 at 1 mM ATP (nM)	: n.a.
