
MSSK1(STK23)

Product code: 04-159

Full-length human MSSK1 [1-567(end) amino acids of accession number NP_055185.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 220
IC50 at 1 mM ATP (nM)	: n.a.

