
MUSK

Product code: 08-153

Human MUSK, catalytic domain [527-869(end) amino acids of accession number NP_005583.1] was expressed as N-terminal GST fusion protein (66 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CSKtide

ATP (μ M) Km app / Bin : 14 / 10
Metal : Mg+Mn
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 2.1
IC50 at 1 mM ATP (nM) : 2.6
