
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 |