

---

---

## SRM(SRMS)

Product code: 08-174

Human SRM, catalytic domain [215-488(end) amino acids of accession number NP\_543013.1] was expressed as N-terminal GST-fusion protein (58kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP ( $\mu$ M) Km app / Bin	: 38 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 290
IC50 at 1 mM ATP (nM)	: 5000