

---

## COT(MAP3K8)

Product code: 07-301

Human COT, catalytic domain [30-397 amino acids of accession number NP\_005195.2] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-COT was purified by using glutathione sepharose chromatography. GST-COT was cleaved by PreScission protease and GST-free COT (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: MAP2K1 Cascade Assay *
ATP ( $\mu$ M)	: 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 120

\* MAP2K1/Erk2/Modified Erktide

---