

COT1/MAP3K8

(Mitogen-activated protein kinase kinase kinase 8, COT proto-oncogene serine/threonine-protein kinase, C-COT, Cancer Osaka thyroid oncogene, M3K8)

- **CAT#:** COT1/MAP3K8
- **Enzyme:** Human COT1/MAP3K8
- **Substrate:** MEK1 (K97R), 1 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Ro 31-8220	17,000
GW 5074	>20,000

- **Scientific Information:** P41279 (<http://www.expasy.org/uniprot/P41279>)

Reference compound IC50 for COT1/MAP3K8

